

17 β -Estradiol ELISA Kit (ADI-900-008)

CELL LYSATE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Melanocytes	Human	M. Fang, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23648481

CULTURE SUPERNATANT

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Vascular smooth muscle cells	Human	A. Bukovski, <i>et al.</i> (2009)	http://www.ncbi.nlm.nih.gov/pubmed/19946214
Embryos	Zebrafish	L. Truong, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27453428

FECES

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Feces	Mixed	T.A. Hanselman, <i>et al.</i> (2004)	http://www.ncbi.nlm.nih.gov/pubmed/15356254
Feces	Octopus	S.E. Larson, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/21192544
Feces	Sea otter	S.E. Larson, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/22753108

PLASMA

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Plasma	Fish	A.K. Sessa, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23651580
Plasma	Fish	C.E. Kidd, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/19607832
Plasma	Human	K. Zhang, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31063987
Plasma	Mouse	P.H. Ear, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30673618
Plasma	Rat	P.H. Ear, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30673618

SERUM

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Serum	Rat	S.K. Kim, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27879453
Serum	Mouse	J.A. Toonen, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27923908
Serum	Mouse	A. Al Mamun, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29692197
Serum	Mouse	J.G. Facciponte, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24642465

TISSUE LYSATE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Ovarian tissue	Mouse	V.K. Maurya, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24628852

WATER

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Aquarium water	Fish	M.R. Kidd, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23168085
Aquarium water	Fish	C.E. Kidd, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/19607832
Seawater	Mixed	S. Heub, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25604269
Manure wastewater	Mixed	T.A. Hanselman, <i>et al.</i> (2004)	http://www.ncbi.nlm.nih.gov/pubmed/15356254
Seawater	Mixed	S. Heub, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25604269

Complete citation can be found on product's webpage

Advance Your Research:
www.enzolifesciences.com

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
 10 Executive Boulevard
 Farmingdale, NY 11735
 Phone: 800.942.0430
 Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
 Industriestrasse 17
 CH-4415 Lausen, Switzerland
 Phone: +41 61 926 8989
 Fax: +41 61 926 8979
info-eu@enzolifesciences.com



Hormone ELISAs: Cited Samples

Progesterone ELISA Kit (ADI-900-011)

CULTURE SUPERNATANT

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Granulosa cells	Cow	T.W. Alemu, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29396041
Endometrial biopsy	Horse	K. Gajos, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25963128
Granulosa cells	Human	Y. Yung, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24147854
Granulosa cells	Human	L. Ophir, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24508664
HEI-193	Human	J.E. Sagers, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29615643
Astrocytes	Mouse	M.A. Mittelman-Smith, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/28384629
LTC-1 cells	Mouse	J.H. Kim, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26739509

FECES

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Feces	Octopus	S.E. Larson, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/21192544

PLASMA

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Plasma	Cow	H. Pothmann, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25670153
Plasma	Fish	L.A. O'Connell, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23899762
Plasma	Fish	L.A. O'Connell, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22166981
Plasma	Fish	C.E. Kidd, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/19607832
Plasma	Horse	J. Kuhl, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26963045
Plasma	Horse	J. Aurich, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30809848
Plasma	Horse	T. Beyer, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30476756
Plasma	Horse	J. Aurich, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25700267
Plasma	Horse	K. Gajos, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25963128
Plasma	Mouse	J. Tanaka, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27335067
Plasma	Mouse	N. Ishii, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31059097
Plasma	Rat	H. Tinwell, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23590819
Plasma	Rat	L.M. Pyter, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23348027
Plasma	Rat	C.M. Contreras, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/28385515
Plasma	Rat	S. Abdulai-Saiku, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/28400143

SERUM

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Serum	Cow	M. Mutinati, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23567219
Serum	Dog	C. Gabriel, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26898415
Serum	Horse	M.J. Siemieniuch, <i>et al.</i> (2016).	https://www.ncbi.nlm.nih.gov/pubmed/27152525
Serum	Mouse	J.G. Facciponte, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24642465
Serum	Mouse	S. Ziegler-Waldkirch, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29855361
Serum	Mouse	T. Nakano, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/29732148
Serum	Pig	A. Dalmau, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25908572
Serum	Rat	E. Estrada-Camarena, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31153863
Serum	Rat	E. Vieyra, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31241201



Hormone ELISAs: Cited Samples

TISSUE LYSATE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Blubber	Dolphin	M.L. Trego, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31110768
Blubber tissue	Whale	L. Pallin, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29942518
Brain	Rat	Z.K. Qiu, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/28740098

WATER

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Aquarium water	Fish	M.R. Kidd, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23168085
Aquarium water	Fish	C.E. Kidd, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/19607832

OTHER SAMPLES

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Follicular fluid	Human	Y. O'Brien, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31217014
Follicular fluid	Pig	P. Pawlak, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30333518
Hair	Brown bear	M. Cattet, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29399362



Hormone ELISAs: Cited Samples

Arg⁸-Vasopressin ELISA Kit (ADI-900-017A)

CULTURE SUPERNATANT

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Neurohypophysis	Rat	T.K. Knott, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18481265

PLASMA

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Plasma	Human	F.K. Idu, <i>et al.</i> (2015)	http://avehjournal.org/index.php/aveh/article/view/10
Plasma	Human	B.A. Corbett, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27540420
Plasma	Human	S.A. Mears, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26805724
Plasma	Human	L.H. Rubin, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27870395
Plasma	Human	O. Oztan, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29309996
Plasma	Human	N. Brondino, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29414026
Plasma	Monkey	R.H. Larke, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27712925
Plasma	Mouse	V. Sauzeau, <i>et al.</i> (2007)	http://www.ncbi.nlm.nih.gov/pubmed/17202406
Plasma	Rat	J.D. Klein, <i>et al.</i> (2006)	http://www.ncbi.nlm.nih.gov/pubmed/16788141
Plasma	Rat	G.E. Haley, <i>et al.</i> (2007)	http://www.ncbi.nlm.nih.gov/pubmed/17522129
Plasma	Rat	G.E. Haley, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18385472
Plasma	Rat	F. Chagnon, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18824919
Plasma	Rat	C. Vega, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/20943859
Plasma	Rat	C. Gray, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23991143
Plasma	Rat	M. Greenwood, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24623760
Plasma	Rat	M.P. Greenwood, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25961839
Plasma	Rat	H.J. Sun, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26746861
Plasma	Rat	A.S. More, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27385784
Plasma	Rat	C.A. Sims, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/29065123
Plasma	Vole	P. Sun, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24561258

SALIVA

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Saliva	Dog	E.L. MacLean, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/28865986
Saliva	Human	O. Weisman, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23246527

SERUM

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Serum	Human	L.H. Rubin, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27870395
Serum	Human	L.H. Rubin, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23465965
Serum	Human	A. Bonavia, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30702630
Serum	Human	L. Dai, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22719898
Serum	Mouse	K. Kaneko, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/21887703
Serum	Rat	B.B. Pinto, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27229490
Serum	Rat	H. Ren, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26517129

TISSUE LYSATE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Brain	Rat	H.F. Zhang, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26265492
Pituitary gland	Rat	M. Greenwood, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24623760
Pituitary gland	Rat	M. Greenwood, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29494864
Pituitary gland	Rat	C.A. Sims, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/29065123

URINE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Urine	Human	P.B. Gray, <i>et al.</i> (2007)	http://www.ncbi.nlm.nih.gov/pubmed/17716675
Urine	Human	C.C. Cohen-Bendahan, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26292177
Urine	Mouse	Y. Zhang, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22622462
Urine	Mouse	Y. Zhang, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26386699
Urine	Mouse	K. Hopp, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25503729
Urine	Mouse	H. Wang, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31033226
Urine	Rat	K. Hopp, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25503729
Urine	Mouse	Y. Zhang, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22622462
Urine	Rat	K. Hopp, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25503729



Hormone ELISAs: Cited Samples

Testosterone ELISA Kit (ADI-900-065)

CELL LYSATE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
LTC-1 cells	Mouse	S.J. Park, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23256993

CEREBROSPINAL FLUID

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Cerebrospinal fluid	Bird	M. Ikeda, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/25342066

CULTURE SUPERNATANT

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Testis explants	Dog	R.G. Lea, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27503122

FECES

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Feces	Gazelle	V.O. Ezenwa, <i>et al.</i> (2011)	http://onlinelibrary.wiley.com/doi/10.1111/j.1365-2435.2011.01919.x/abstract
Feces	Octopus	S.E. Larson, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/21192544

PLASMA

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Plasma	Bird	M.S. DeVries, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/21986087
Plasma	Bird	S. Davies, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26163581
Plasma	Bird	S. Davies, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25985895
Plasma	Bird	S.E. Lynn, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25849310
Plasma	Bird	C.M. Bauer, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26718082
Plasma	Bird	P. Deviche, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27311790
Plasma	Bird	H.E. Watts, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26836771
Plasma	Bird	K.B. Needham, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/28017731
Plasma	Bird	S. Davies, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29166572
Plasma	Bird	I.I. Levin, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30341385
Plasma	Bird	A.R. Robart, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30225078
Plasma	Bird	J.L. Graham, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31236557
Plasma	Fish	A.K. Greenwood, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18628117
Plasma	Fish	C.E. Kidd, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/19607832
Plasma	Fish	L.A. O'Connell, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22166981
Plasma	Fish	A.K. Sessa, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23651580
Plasma	Horse	A. Lemasson, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25714814
Plasma	Lizard	R.J. Seddon, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27137079
Plasma	Mouse	P. Manna, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18926901
Plasma	Mouse	T. Homma, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27888021
Plasma	Mouse	C. Fu, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29728588
Plasma	Mouse	H.H. Shawki, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29324782
Plasma	Rabbit	A.M. Traish, <i>et al.</i> (2003)	http://www.ncbi.nlm.nih.gov/pubmed/12721214
Plasma	Rat	H. Tinwell, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23590819



Hormone ELISAs: Cited Samples

Plasma	Rat	A.P. Borrow, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23419048
Plasma	Rat	M. Choi, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23548559
Plasma	Rat	V. Diwan, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/25209863
Plasma	Rat	N. Yamaguchi, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27981176
Plasma	Rat	J.H. Choi, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27550492
Plasma	Rat	V. Brouard, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27174447
Plasma	Rat	S. Abdulai-Saiku, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/28400143
Plasma	Rat	A. Asimes, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30841593
Plasma	Sea turtle	M.P. Jensen, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29316410

SERUM

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Serum	Bat	A.C.F. Souza, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30332444
Serum	Fish	A.K. Singh, <i>et al.</i> (2015)	http://www.omicsonline.com/open-access/a-cyp-based-sex-determination-and-monosex-production-in-aquaculture-species-oreochromis-niloticus-l-and-a-cyprinid-cyprinus-carpio-l2150-3508.10000112.php?aid=38985
Serum	Hamster	J.C. Wommack, <i>et al.</i> (2004)	http://www.ncbi.nlm.nih.gov/pubmed/15344915
Serum	Hamster	G.E. Demas, <i>et al.</i> (2004)	http://www.ncbi.nlm.nih.gov/pubmed/15555500
Serum	Hamster	T.J. Greives, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/19094081
Serum	Hamster	K.L.P. Long, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30297514
Serum	Mole-rat	D.E. Peragine, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27018426
Serum	Mouse	C. Chang, <i>et al.</i> (2004)	http://www.ncbi.nlm.nih.gov/pubmed/15107499
Serum	Mouse	A. de Peyster, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18824092
Serum	Mouse	R.N. Thompson, <i>et al.</i> (2015)	http://www.omicsgroup.org/journals/prenatal-testosterone-exposure-influences-neuronal-sensitivity-to-pheromones-in-female-mice-2168-9652-1000170.php?aid=59811
Serum	Mouse	M.S. Latsko, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27108196
Serum	Mouse	J.A. Toonen, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27923908
Serum	Mouse	A. Al Mamun, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29692197
Serum	Mouse	Y. Hu, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/28982275
Serum	Mouse	H.J. Lee, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29689290
Serum	Mouse	J.S. Templin, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31047887
Serum	Mouse	B. Weger, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30344015
Serum	Rabbit	A. Panossian, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/21901061
Serum	Rat	D.L. Cyrenne, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/21920363
Serum	Rat	J.H. Choi, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27550492
Serum	Rat	J.F. Lynch 3rd, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27368147
Serum	Rat	J.M. Moon, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/28830494
Serum	Rat	C.T. Ribeiro, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29384140
Serum	Rat	C.T. Ribeiro, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30600770
Serum	Rat	V. Migliaccio, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31083466



Hormone ELISAs: Cited Samples

TISSUE LYSATE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Antler	Velvet	S.H. Tseng, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22951396
Blubber	Whale	K.A. Cates, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30904390
Hair	Brown bear	M. Cattet, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29399362
Hair	Rat	L. Arnon, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27125699
Mixed	Snail	N.E. Omran, <i>et al.</i> (2016)	http://www.ncbi.nlm.nih.gov/pubmed/24215068
Mixed	Snail	A.M. Ibrahim, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30040958
Testis	Mouse	K. Ko, <i>et al.</i> (2002)	http://www.ncbi.nlm.nih.gov/pubmed/12441367
Testis	Rat	M. Clarke, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24499511

WATER

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Aquarium water	Fish	M.R. Kidd, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23168085
Aquarium water	Fish	C.E. Kidd, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/19607832

OTHER SAMPLES

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Hemolymph	Snail	A.M. Ibrahim, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30684178
Sperm	Rat	C.T. Ribeiro, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29384140



Hormone ELISAs: Cited Samples

Cortisol ELISA Kit (ADI-900-071)

AMNIOTIC FLUID

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Amniotic fluid	Pig	I. Trus, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30058440
Amniotic fluid	Rat	Y. Jang, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29183725

CULTURE SUPERNATANT

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Caco-2 cells	Human	Y. Tomabechei, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/28744834
H295R cells	Human	T. Escajadillo, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26992564
H295R cells	Human	N.C. Lucki, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/21864647
Liver and fat pads	Mouse	H. Ho, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26519792
Liver and fat pads	Mouse	S.B. Park, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26515507

FECES

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Feces	Gazelle	V.O. Ezenwa, <i>et al.</i> (2011)	http://onlinelibrary.wiley.com/doi/10.1111/j.1365-2435.2011.01919.x/abstract
Feces	Monkey	C. Cinque, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27904967
Feces	Mouse	A. Running (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26170887

MILK

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Milk	Human	L. O'Rourke, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29686862

PLASMA

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Plasma	Cow	M.E. Olson, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27295955
Plasma	Cow	B. Earley, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26520781
Plasma	Cow	A.K. Kelly, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/28804620
Plasma	Fish	J.H. Kim, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26991846
Plasma	Fish	B.C. Mommer, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23628383
Plasma	Fish	C.A. Wei, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/28734797
Plasma	Fish	J.H. Kim, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27689810
Plasma	Fish	P.J. Tapia, <i>et al.</i> (2012)	http://www.tandfonline.com/doi/abs/10.1080/15222055.2012.676020?journalCode=unaj20
Plasma	Fish	A.K. Greenwood, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18628117
Plasma	Horse	A. Lemasson, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25714814
Plasma	Human	P.W. Watt, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27157340
Plasma	Human	E. Skornyakov, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30089765
Plasma	Human	L.A. Ray, <i>et al.</i> (2009)	http://www.ncbi.nlm.nih.gov/pubmed/18824022
Plasma	Pig	J.L. Rault, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26479373
Plasma	Pig	D.L. Turpin, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27792138
Plasma	Pig	D.L. Turpin, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26999224
Plasma	Pig	I. Trus, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30058440
Plasma	Rat	Y. Jang, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29183725

TISSUE LYSATE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Tissue	Fish	G.D. King, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26780133
Epididymal fat pads	Mouse	H. Oh, <i>et al.</i> (2015)	https://www.ncbi.nlm.nih.gov/pubmed/26519792
Lamina	Whale	S.J. Trumble, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30389921
Liver	Mouse	H. Oh, <i>et al.</i> (2015)	https://www.ncbi.nlm.nih.gov/pubmed/26519792
Whole body tissue	Fish larvae	G. Luzeena Raja, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30395803

SALIVA

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Saliva	Horse	J. Aurich, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25700267
Saliva	Human	L. Starnino, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27788238
Saliva	Human	A.P. Allen, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27801892
Saliva	Human	E.A. Nelson, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18789315
Saliva	Pig	M. Zupan, <i>et al.</i> (2017)	http://www.scielo.br/scielo.php?script=sci_arttext&pid=S1516-35982017000100033

SERUM

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Serum	Fish	N.A. Khan, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30847628
Serum	Hamster	E.D. Carlton, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26542473
Serum	Human	A.W. Carrico, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29266420
Serum	Hamster	J.C. Wommack, <i>et al.</i> (2004)	http://www.ncbi.nlm.nih.gov/pubmed/15344915
Serum	Hamster	G.E. Demas, <i>et al.</i> (2004)	http://www.ncbi.nlm.nih.gov/pubmed/15555500
Serum	Hamster	J.C. Wommack, <i>et al.</i> (2003)	http://www.ncbi.nlm.nih.gov/pubmed/14637229
Serum	Mouse	C. Raja Gabaglia, <i>et al.</i> (2007)	http://www.ncbi.nlm.nih.gov/pubmed/17332360
Serum	Mouse	N.R. Han, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29286166
Serum	Rabbit	W.P. Cawthorn, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26696121
Serum	Rabbit	A. Panossian, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/21901061

URINE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Urine	Human	H. Fukuda, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/19146551

WATER

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Aquarium water	Fish	L.J. Henderson, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/29018617



Hormone ELISAs: Cited Samples

Corticosterone ELISA Kit (ADI-900-097)

AMNIOTIC FLUID

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Amniotic fluid	Mouse	X. Xiang, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27974238

CULTURE SUPERNATANT

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Melanocytes	Human	A. Slominski, <i>et al.</i> (2005)	http://www.ncbi.nlm.nih.gov/pubmed/15572653
Y-1 adrenal cells	Mouse	M.S. Charles, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25816736

FECES

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Feces	Lynx	K.V. Fanson, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/21717144
Feces	Mouse	R.I. Monarca, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26143188
Feces	Mouse	H. Siemen, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/22016787
Feces	Rat	P.V. Turner, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22330864
Feces	Rat	A. Green, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29183765
Feces	Rat	J.D. Vega-Torres, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29428401
Feces	Rat	E.D.K. Sullivan, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30708127

PLASMA

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Plasma	Alligator	J.W. Finger, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30976886
Plasma	Alligator	M.T. Hamilton, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30031024
Plasma	Bird	O.P. Love, <i>et al.</i> (2005)	http://www.ncbi.nlm.nih.gov/pubmed/16475090
Plasma	Bird	T. Tachibana, <i>et al.</i> (2007)	http://www.ncbi.nlm.nih.gov/pubmed/17321768
Plasma	Bird	G.S. Archer, <i>et al.</i> (2015)	http://scialert.net/abstract/?doi=ijps.2015.293.299
Plasma	Bird	R.J. Steenweg, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25913259
Plasma	Bird	V. Marasco, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26093051
Plasma	Bird	W.A. Buttemer, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25701624
Plasma	Bird	S.E. Lynn, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25849310
Plasma	Bird	J.C. Huth, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26201348
Plasma	Bird	C. Eikenaar, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26163918
Plasma	Bird	P. Deviche, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27311790
Plasma	Bird	H.L. Hennin, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26925215
Plasma	Bird	O.L. Crino, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26828818
Plasma	Bird	Y. Aharon-Rotman, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/26699204
Plasma	Bird	S. Gao, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29530549
Plasma	Bird	O.L. Crino, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29197553
Plasma	Chicken	K. Furukawa, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26883687
Plasma	Bird	S. Davies, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29166572
Plasma	Bird	C. Eikenaar, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29958878
Plasma	Bird	S. Guindre-Parker, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30190315
Plasma	Bird	I.I. Levin, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30341385
Plasma	Bird	B. Jimeno, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30158537
Plasma	Bird	B. Jimeno, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30322980



Hormone ELISAs: Cited Samples

Plasma	Bird	J.R. Nelson, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30287758
Plasma	Bird	A.R. Robart, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30225078
Plasma	Bird	L.N. Cooper, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30517055
Plasma	Bird	K. Sarpong, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30392885
Plasma	Chicken	M. Ericsson, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27105229
Plasma	Chicken	F. Gualtieri, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31073135
Plasma	Chicken	L. Hedlund, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30787406
Plasma	Fish	L.A. O'Connell, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23899762
Plasma	Human	Q.R. Xia, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29427582
Plasma	Lizard	R.J. Seddon, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27137079
Plasma	Mouse	Z. Zhai, <i>et al.</i> (2009)	http://www.ncbi.nlm.nih.gov/pubmed/19303756
Plasma	Mouse	P.A. Volden, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23780289
Plasma	Mouse	Z. Liu, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26133658
Plasma	Mouse	X. Liu, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25959659
Plasma	Mouse	L.C. Lin, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25600109
Plasma	Mouse	C. Yamada, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25732068
Plasma	Mouse	E. Lara-Padilla, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26198915
Plasma	Mouse	F. Reichmann, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26066467
Plasma	Mouse	H.M. Savignac, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25794930
Plasma	Mouse	S.N. Batchu, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25999402
Plasma	Mouse	C.Y. Tsai, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26415720
Plasma	Mouse	K. Ramirez, <i>et al.</i> (2016)	http://www.ncbi.nlm.nih.gov/pubmed/26342944
Plasma	Mouse	T.H. Tran, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26850477
Plasma	Mouse	N. Soualeh, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27702682
Plasma	Mouse	G.R. Rompala, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27286933
Plasma	Mouse	S. Chang, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27701413
Plasma	Mouse	R. Chavan, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26838474
Plasma	Mouse	J.M. Castellano, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27364522
Plasma	Mouse	A. Cabral, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26874559
Plasma	Mouse	K. Ramirez, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26342944
Plasma	Mouse	S. Arase, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27500935
Plasma	Mouse	R. Mayerhofer, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27751870
Plasma	Mouse	D. Caballero-Hernandez, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/29284778
Plasma	Mouse	Y.A. Lee, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29472615
Plasma	Mouse	L.M. Pyter, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29229147
Plasma	Mouse	K.H. Seo, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29928889
Plasma	Mouse	K.M. Corder, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30024942
Plasma	Mouse	A. Niraula, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29382712
Plasma	Mouse	C.D. Staton, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30240784
Plasma	Mouse	H. Zhao, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30003515
Plasma	Mouse	S. Abdurakhmanova, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31216095
Plasma	Mouse	A.V. Aubry, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30542090
Plasma	Mouse	X. Gorny, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30797047
Plasma	Mouse	H. Li, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31241809



Hormone ELISAs: Cited Samples

Plasma	Mouse	K.A. Sullivan, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31108169
Plasma	Rat	S. Ellis, <i>et al.</i> (2005)	http://www.ncbi.nlm.nih.gov/pubmed/15972802
Plasma	Rat	L. Desbonnet, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18456279
Plasma	Rat	S.M. McBride, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18614766
Plasma	Rat	Q.Y. Meng, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/21664419
Plasma	Rat	C.M. O'Mahony, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/21110995
Plasma	Rat	B. Ray, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/21477639
Plasma	Rat	M.S. Charles, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22779090
Plasma	Rat	N. Lajud, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/21862224
Plasma	Rat	M.S. Khalil (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26379312
Plasma	Rat	P.K. Wang, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26342649
Plasma	Rat	N.C. Donner, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26454419
Plasma	Rat	M.R. Favre, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26089770
Plasma	Rat	C.M. Sirieix, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26059034
Plasma	Rat	G.S. Padzys, <i>et al.</i> (2015)	http://peertechz.blogspot.ch/2015/10/impact-of-early-nasal-obstruction-in.html
Plasma	Rat	J. Bourgin, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25823695
Plasma	Rat	C.S. Harrell, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26356038
Plasma	Rat	K.E. Boschen, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25805052
Plasma	Rat	Y. Ruvalcaba-Delgadillo, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26168952
Plasma	Rat	H. Er, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25596037
Plasma	Rat	G. Leo, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25654993
Plasma	Rat	M. Bülbül, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27829122
Plasma	Rat	A. Gano, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27208497
Plasma	Rat	G.G. Page, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27981183
Plasma	Rat	A.B. Jones, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27247143
Plasma	Rat	H.J. Lee, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26903707
Plasma	Rat	E.A. Karanges, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26876759
Plasma	Rat	M. Bekhbat, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26946276
Plasma	Rat	A.L. Feldhaus, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27906551
Plasma	Rat	J.H. Choi, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27550492
Plasma	Rat	W.E. Patarroyo, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27913253
Plasma	Rat	A. Torcaso, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27930953
Plasma	Rat	W. Wang, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/28296942
Plasma	Rat	C.M. Contreras, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/28385515
Plasma	Rat	M. Bekhbat, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29730284
Plasma	Rat	A.C. Kentner, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30121011
Plasma	Rat	V. Satta, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29456490
Plasma	Rat	M.H. Ali, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30954669
Plasma	Rat	M. Bekhbat, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30550932
Plasma	Rat	A.S.M. Hasan Mahmood, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30660767
Plasma	Rat	D. Huzard, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30448728
Plasma	Rat	S.K. Mandal, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30481553
Plasma	Toad	S. Gardner, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30030104
Plasma	Turtle	J.M. West, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29548757



Hormone ELISAs: Cited Samples

Plasma	Vole	M.A. Scotti, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25639952
Plasma	Vole	J.R. Yee, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26836772
Plasma	Vole	N. McNeal, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/28008793
Plasma	Vole	M. Soto, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30769014

SERUM

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Serum	Chicken	P.A. Kolodziejski, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29416857
Serum	Chicken	S. Gurung, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29324639
Serum	Mouse	J. Penney, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27789393
Serum	Mouse	O.K. Sial, <i>et al.</i> (2016)	http://www.ncbi.nlm.nih.gov/pubmed/26545443
Serum	Mouse	K. Saito, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26988598
Serum	Mouse	C.F. Li, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27018164
Serum	Mouse	M.S. Latsko, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27108196
Serum	Mouse	K.S. Abelson, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27815456
Serum	Mouse	H. Liu, <i>et al.</i> (2015)	http://www.scirp.org/journal/PaperInformation.aspx?Paper-ID=60373
Serum	Mouse	J. Burkus, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25985793
Serum	Mouse	S. Shushimita, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26616751
Serum	Mouse	Z.Z. Wang, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26161529
Serum	Mouse	Y. Zhang, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25603858
Serum	Mouse	Y. Mishima, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25754523
Serum	Mouse	S. Yokota, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26291742
Serum	Mouse	L. Desbonnet, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25866195
Serum	Mouse	M. Ghilan, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25731748
Serum	Mouse	P. Wong, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25964750
Serum	Mouse	N.L. Baganz, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26529424
Serum	Mouse	H. Siemen, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/22016787
Serum	Mouse	W.L. Thompson, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18706086
Serum	Mouse	A.P. Fernandez, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18723674
Serum	Mouse	M. Aghajani, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/29145213
Serum	Mouse	J.E. Choi, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27988398
Serum	Mouse	G. Caratti, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30179226
Serum	Mouse	K.S. Collison, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29561882
Serum	Mouse	S. Moreno, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30127743
Serum	Mouse	M.A. Quinn, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29514097
Serum	Mouse	H. Woo, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29970188
Serum	Mouse	H. Baazim, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31110314
Serum	Mouse	N. Dos Santos, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31040362
Serum	Mouse	L.M. Ince, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/29965797
Serum	Mouse	I.H.N. Luijten, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31067456
Serum	Mouse	J. Shah, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30691851
Serum	Rabbit	W.P. Cawthorn, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26696121
Serum	Rat	L.R. Jacinto, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27940019
Serum	Rat	L.S. Novaes, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27815155



Hormone ELISAs: Cited Samples

Serum	Rat	L.M. Zhang, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27845056
Serum	Rat	L.L. Williamson, <i>et al.</i> (2016)	http://www.ncbi.nlm.nih.gov/pubmed/26162711
Serum	Rat	A. Gano, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27208497
Serum	Rat	A.S. Koe, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26836417
Serum	Rat	Y.E. Lin, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27109341
Serum	Rat	D. Bonini, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26966584
Serum	Rat	R. Schneider, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25771148
Serum	Rat	P. Licznarski, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26506154
Serum	Rat	M.H. Nowland, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/21535924
Serum	Rat	A.P. Johnston, <i>et al.</i> (2007)	http://www.ncbi.nlm.nih.gov/pubmed/17092995
Serum	Rat	L.K. Fonken, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29524458
Serum	Rat	S. Sun, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30110343
Serum	Rat	P. Chakraborty, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31193585
Serum	Rat	M. Navarro-Meza, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31060976
Serum	Rat	Q. Pan, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30699999
Serum	Rat	N. Zhao, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31201489

TISSUE LYSATE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Claw	Chameleon	D. Matas, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26993343
Hippocampus	Mouse	J.S. Bae, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27792185
Amygdala	Rat	M.G. Frank, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31078691
Adrenal gland	Mouse	N.R. Mahapatra, <i>et al.</i> (2005)	http://www.ncbi.nlm.nih.gov/pubmed/16007257
Hair	Rat	L. Arnon, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27125699
Hippocampus	Rat	J.L. Sobesky, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27595136
Hippocampus	Rat	B. Bray, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27208490
Hippocampus	Rat	M.G. Frank, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/31078691
Scute	Alligator	M.T. Hamilton, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30031024

URINE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Urine	Mouse	E.K. Hutchinson, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22330863
Urine	Mouse	A.E. Fitchett, <i>et al.</i> (2005)	http://www.ncbi.nlm.nih.gov/pubmed/16098684

WHOLE BLOOD

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Blood	Mouse	P. Mancuso, <i>et al.</i> (2006)	http://www.ncbi.nlm.nih.gov/pubmed/16210671



Hormone ELISAs: Cited Samples

Oxytocin ELISA Kit (ADI-900-153A)

CEREBROSPINAL FLUID

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Cerebrospinal fluid	Human	C.L. Clark, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/23507187
Cerebrospinal fluid	Monkey	S.M. Freeman, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26826355
Cerebrospinal fluid	Rat	N.S. Couto-Pereira, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27717870
Cerebrospinal fluid	Mouse	A. Oppong-Damoah, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30593782
Cerebrospinal fluid	Rat	G. Martinez-Lorenzana, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18823708

PLASMA

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Plasma	Dog	M. Petersson, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/29081760
Plasma	Human	L.H. Rubin, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27870395
Plasma	Human	V.M. Husarova, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27081377
Plasma	Human	H. Kosaka, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27552585
Plasma	Human	B.A. Corbett, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27540420
Plasma	Human	R.T. Emeny, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25827957
Plasma	Human	S.H. Lin, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25871910
Plasma	Human	D.A. Muin, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26151620
Plasma	Human	M. Eidelman-Rothman, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26175673
Plasma	Human	P.Y. Lin, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23460821
Plasma	Human	A. Szeto, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/21636661
Plasma	Human	R. Feldman, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/20558167
Plasma	Human	J. Holt-Lunstad, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18842740
Plasma	Human	O. Oztan, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29309996
Plasma	Human	N. Brondino, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29414026
Plasma	Human	M. Petersson, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/29081760
Plasma	Human	N. Kampka, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30539300
Plasma	Monkey	R.H. Larke, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27712925
Plasma	Monkey	S.M. Freeman, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26826355
Plasma	Mouse	O. Lopatina, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/21501260
Plasma	Mouse	N. Scott, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26375004
Plasma	Mouse	P.H. Ear, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30673618
Plasma	Mouse	A. Oppong-Damoah, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30593782
Plasma	Pig	M. Norrby, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/20626675
Plasma	Pig	M. Marcet Rius, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29353556
Plasma	Rabbit	A. Szeto, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/22998949
Plasma	Rat	M. Bülbül, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27829122
Plasma	Rat	A.S. Pereira, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26297864
Plasma	Rat	L.M. Carini, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/24005186
Plasma	Rat	C. Vega, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/20943859
Plasma	Rat	G.E. Haley, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18385472
Plasma	Rat	G. Martinez-Lorenzana, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18823708
Plasma	Rat	G.E. Haley, <i>et al.</i> (2007)	http://www.ncbi.nlm.nih.gov/pubmed/17522129
Plasma	Rat	P.H. Ear, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30673618
Plasma	Rat	A. Holubova, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30984017



Hormone ELISAs: Cited Samples

Plasma	Seal	K.J. Robinson, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26246656
Plasma	Vole	J.R. Yee, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26836772
Plasma	Vole	P. Sun, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24561258

SALIVA

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Saliva	Dog	E.L. MacLean, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/28865986
Saliva	Human	S.M. Ormsby, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26887958
Saliva	Human	G.A. Behr, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27989492
Saliva	Human	A.V. Jaeggi, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25788487
Saliva	Human	S. Tsuji, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25954210
Saliva	Human	E.R. Lebowitz, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26716876
Saliva	Human	R. Feldman, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/20558167
Saliva	Human	J. Holt-Lunstad, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18842740
Saliva	Human	K. Takahata, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29447299

SERUM

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Serum	Dog	F.K. Hollinshead, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/20171720
Serum	Human	T. Sasaki, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27536785
Serum	Human	L. Dai, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22719898
Serum	Human	L.H. Rubin, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/23465965
Serum	Human	V.B. Morhenn, <i>et al.</i> (2008)	http://www.ehonline.org/article/S1090-5138(08)00048-2/abstract
Serum	Human	G.A. Behr, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/27989492
Serum	Monkey	V. Michopoulos, <i>et al.</i> (2011)	http://www.ncbi.nlm.nih.gov/pubmed/21316367
Serum	Mouse	J. Roizen, <i>et al.</i> (2007)	http://www.ncbi.nlm.nih.gov/pubmed/17895964

TISSUE LYSATE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Brain	Rat	H.F. Zhang, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26265492
Olfactory bulb	Rat	A.C. Meidahl, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29266199
Pons	Rat	A.C. Meidahl, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29266199
Spinal cord	Rat	A.C. Meidahl, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29266199
Lumbar spinal cord	Rat	G. Martinez-Lorenzana, <i>et al.</i> (2008)	http://www.ncbi.nlm.nih.gov/pubmed/18823708
Trigeminal ganglia	Rat	A.C. Meidahl, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29266199



Hormone ELISAs: Cited Samples

URINE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Urine	Human	R. Gregory, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25562711
Urine	Human	C.C. Cohen-Bendahan, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26292177
Urine	Human	K.G. Salvante, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22121074
Urine	Human	J. Bick, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/20953313
Urine	Human	P.B. Gray, <i>et al.</i> (2007)	http://www.ncbi.nlm.nih.gov/pubmed/17716675
Urine	Human	S.B. Algoe, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/28968183
Urine	Human	Y. Eisenberg, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29300770
Urine	Monkey	L.R. Moscovice, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22986337
Urine	Monkey	C. Finkenwirth, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26232089
Urine	Monkey	C. Finkenwirth, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26836769
Urine	Tamarin	C.T. Snowdon, <i>et al.</i> (2010)	http://www.ncbi.nlm.nih.gov/pubmed/20600045



Hormone ELISAs: Cited Samples

Aldosterone ELISA Kit (ADI-900-173)

PLASMA

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Plasma	Dog	X. Wang, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25349225
Plasma	Human	S.A. Mears, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/26805724
Plasma	Human	V. Conway, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24206823
Plasma	Monkey	A.L. Feldhaus, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27906551
Plasma	Mouse	C.K. Hathaway, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/25902541
Plasma	Mouse	Y.N. Qiao, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24951589
Plasma	Rat	A.S. More, <i>et al.</i> (2016)	https://www.ncbi.nlm.nih.gov/pubmed/27385784
Plasma	Rat	I. Dhar, <i>et al.</i> (2014)	http://www.ncbi.nlm.nih.gov/pubmed/24436324
Plasma	Rat	A. Mansuri, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/29051307
Plasma	Rat	A.M.A. de Souza, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29985423
Plasma	Rat	G. Zhou, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29848510

SERUM

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Serum	Mouse	J. Mittag, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23257356
Serum	Mouse	H. Li, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/29499204
Serum	Rat	S. Sakata, <i>et al.</i> (2015)	http://www.scirp.org/journal/PaperInformation.aspx?paperID=58692
Serum	Rat	S.K. Singh, <i>et al.</i> (2012)	http://www.ncbi.nlm.nih.gov/pubmed/22322970
Serum	Rat	S. Nagao, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30870430

URINE

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Urine	Mouse	Y. Zhang, <i>et al.</i> (2015)	http://www.ncbi.nlm.nih.gov/pubmed/26386699
Urine	Rat	T. Katayama, <i>et al.</i> (2013)	http://www.ncbi.nlm.nih.gov/pubmed/23974905
Urine	Rat	D.V. Ilatovskaya, <i>et al.</i> (2019)	https://www.ncbi.nlm.nih.gov/pubmed/30745171

OTHER SAMPLES

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Kidney	Mouse	Z. Li, <i>et al.</i> (2018)	https://www.ncbi.nlm.nih.gov/pubmed/30623139
Vitreous humor	Rat	M. Kobayashi, <i>et al.</i> (2017)	https://www.ncbi.nlm.nih.gov/pubmed/28253402