



Wildlife Endocrinology: Cited Sample Types

Aldosterone ELISA Kit (ADI-900-173)

SPECIES	SAMPLE ORIGIN	REFERENCES
Dog	Plasma	Wang et al., 2015
Human	Plasma	Mears et al., 2016 ; Conway et al., 2014
Mouse	Urine	Zhang et al., 2015
Mouse	Plasma	Hathaway et al., 2015 ; Qiao et al., 2014
Mouse	Serum	Mittag et al., 2013
Rat	Serum	Sakata et al., 2015
Rat	Plasma	Dhar et al., 2014
Rat	Urine	Katatayama et al., 2013

Angiotensin II ELISA Kit (ADI-900-204)

SPECIES	SAMPLE ORIGIN	REFERENCES
Human	Tissue Lysate	Holappa et al., 2015
Mouse	Tissue Lysate	Werner et al., 2016
Mouse	Plasma	Chen et al., 2014 ; Tsukuda et al., 2012
Mouse	Serum	Mittag et al., 2013
Pig	Plasma	Jakob et al., 2012
Rat	Tissue Lysate	Xu et al., 2016
Rat	Plasma	Waghe et al., 2015 ; Park et al., 2015

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

SPECIES	SAMPLE ORIGIN	REFERENCES
Human	Serum	Mears et al., 2016 ; Rubin et al., 2013 ; Dai et al., 2012
Human	Urine	Cohen-Bendahan et al., 2015 ; Gray et al., 2007
Human	Plasma	Idu et al., 2015
Human	Saliva	Weisman et al., 2013
Mouse	Urine	Zhang et al., 2015 ; Hopp et al., 2015 ; Zhang et al., 2012
Mouse	Serum	Kaneko et al., 2011
Mouse	Plasma	Sauzeau et al., 2007
Rat	Plasma	Pinto et al., 2016 ; Sun et al., 2016 ; Greenwood et al., 2015 ; Greenwood et al., 2014 ; Gray et al., 2013 ; Vega et al., 2010 ; Chagnon et al., 2008 ; Haley and Flynn, 2008 ; Haley and Flynn, 2007 ; Klein et al., 2006
Rat	Urine	Hopp et al., 2015
Rat	Tissue Lysate	Zhang et al., 2015 ; Greenwood et al., 2014
Rat	Serum	Ren et al., 2015
Rat	Culture Supernatant	Knott et al., 2008
Vole	Plasma	Sun et al., 2014

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

Corticosterone ELISA Kit (ADI-900-097)

SPECIES	SAMPLE ORIGIN	REFERENCES
Bird	Plasma	Crino et al., 2016 ; Hennin et al., 2016 ; Deviche et al., 2016 ; Lynn et al., 2015 ; Aharon-Rotman et al., 2015 ; Eikenaar et al., 2015 ; Buttemer et al., 2015 ; Deviche et al., 2014 ; Mark and Rubenstein, 2013 ; Love et al., 2005
Chameleon	Tissue Lysate	Matas et al., 2016
Chicken	Plasma	Ericsson et al., 2016 ; Furukawa et al., 2016 ; Huth and Archer, 2015 ; Archer, 2015
Duck	Plasma	Steenweg et al., 2015
Fish	Plasma	O'Connell et al., 2013
Human	Culture Supernatant	Slominski et al., 2005
Lizard	Plasma	Seddon and Hews, 2016
Lynx	Feces	Fanson et al., 2012
Mouse	Serum	Latsko et al., 2016 ; Li et al., 2016 ; Cawthorn et al., 2016 ; Saito et al., 2016 ; Shushimita et al., 2016 ; Wang et al., 2015 ; Ghilan et al., 2015 ; Desbonnet et al., 2015 ; Yokota et al., 2015 ; Mishima et al., 2015 ; Zhang et al., 2015 ; Wong et al., 2015 ; Burkus et al., 2015 ; Liu et al., 2015 ; Sial et al., 2015
Mouse	Plasma	Ramirez et al., 2016 ; Cabral et al., 2016 ; Chavan et al., 2016 ; Rompala et al., 2016 ; Tran et al., 2016 ; Baganz et al., 2015 ; Tsai et al., 2015 ; Batchu et al., 2015 ; Savignac et al., 2015 ; Reichmann et al., 2015 ; Lara-Padilla et al., 2015 ; Yamada et al., 2015 ; Lin and Sibille, 2015 ; Lui et al., 2015 ; Vega-Rivera et al., 2015 ; Lui et al., 2015 ; Zhai et al., 2009 ; Fernandez et al., 2008
Mouse	Feces	Siemen et al., 2011
Mouse	Tissue Lysate	Mahapatra et al., 2005
Mouse	Urine	Hutchinson et al., 2012 ; Fitchett et al., 2005
Rat	Serum	Bonini et al., 2016 ; Lin et al., 2016 ; Williamson et al., 2016 ; Koe et al., 2016 ; Licznernski et al., 2015 ; Schneider et al., 2015 ; Wang et al., 2015 ; Nowland et al., 2011 ; Johnston et al., 2007
Rat	Plasma	Bray et al., 2016 ; Bekhbat et al., 2016 ; Karanges et al., 2016 ; Lee et al., 2016 ; Leo et al., 2015 ; Er et al., 2015 ; Ruvalcaba-Delgado et al., 2015 ; Boschen et al., 2015 ; Harrell et al., 2015 ; Bourgin et al., 2015 ; Padzys et al., 2015 ; Sirieix et al., 2015 ; Favre et al., 2015 ; Donner et al., 2015 ; Khalil, 2015 ; Lajud et al., 2012 ; Charles et al., 2012 ; Aydin et al., 2011 ; O'Mahony et al., 2011 ; Meng et al., 2011 ; McBride et al., 2008 ; Desbonnet et al., 2008 ; Ellis et al., 2005
Rat	Hair	Arnon et al., 2016
Rat	Feces	Turner et al., 2012
Vole	Plasma	Yee et al., 2016 ; Scotti et al., 2015

Cortisol ELISA Kit (ADI-900-071)

SPECIES	SAMPLE ORIGIN	REFERENCES
Cow	Plasma	Olson et al., 2016 ; Earley et al., 2015
Fish	Tissue Lysate	King et al., 2016
Fish	Plasma	Kim and Kang, 2016 ; Mommer and Bell, 2013 ; Tapia et al., 2012 ; Greenwood et al., 2008
Gazelle	Feces	Ezenwa et al., 2012
Hamster	Serum	Carlton and Demas, 2015 ; Demas et al., 2004 ; Wommack and Delville, 2003
Hamster	Plasma	Wommack et al., 2004
Horse	Plasma	Lemasson et al., 2015
Human	Plasma	Watt et al., 2016
Human	Culture Supernatant	Lucki et al., 2012
Human	Saliva	Nelson et al., 2008
Human	Urine	Fukuda et al., 2008
Mouse	Serum	Cawthorn et al., 2016 ; Gabaglia et al., 2007
Mouse	Tissue Lysate	Oh et al., 2015
Mouse	Feces	Running, 2015
Pig	Plasma	Turpin et al., 2016 ; Rault et al., 2015

Direct 8-ISO-PGF_{2α} ELISA Kit (ADI-900-091)

SPECIES	SAMPLE ORIGIN	REFERENCES
Dog	Plasma	Hinchcliff et al., 2000
Hamster	Culture Supernatant	Patel and Bhat, 2004
Hamster	Tissue Lysate	Bhat et al., 2003
Human	Tissue Lysate	Johnston et al., 2016
Human	Plasma	Duffy et al., 2015 ; Kumar et al., 2015 ; Li et al., 2014 ; Jeanes et al., 2004
Human	Serum	Puthuchery and Nathan, 2008
Human	Urine	Huang et al., 2002
Human	Culture Supernatant	Patel and Bhat, 2004
Mouse	Plasma	Hoffman-Goetz and Quadrilatero, 2003
Rat	Tissue Lysate	Fang and Lin, 2008

DHEA ELISA Kit (ADI-900-093)

SPECIES	SAMPLE ORIGIN	REFERENCES
Cow	Plasma	O'Driscoll et al., 2015
Human	Serum	Sato et al., 2016
Human	Culture Supernatant	Escajadillo et al., 2016 ; Lucki et al., 2012

17β-Estradiol ELISA Kit (ADI-900-008)

SPECIES	SAMPLE ORIGIN	REFERENCES
Fish	Plasma	Sessa et al., 2013
Human	Culture Supernatant	Fang et al., 2013 ; Bukovsky, 2009
Mouse	Tissue Lysate	Maurya et al., 2014
Mouse	Serum	Facciponte et al., 2014
Octopus	Feces	Larson and Anderson, 2010
Sea otter	Feces	Larson et al., 2013
	Water	Heub et al., 2015 ; Kidd et al., 2013 ; Kidd et al., 2010 ; Hanselman et al., 2004

17β-Estradiol High Sensitivity ELISA Kit (ADI-900-174)

SPECIES	SAMPLE ORIGIN	REFERENCES
Cow	Serum	Mutinati et al., 2013
Fish	Serum	Singh and Srivastava, 2015
Gerbil	Tissue Lysate	Zanatelli et al., 2014
Goat	Plasma	Castagnino et al., 2015
Human	Plasma	Zolin et al., 2015
Mouse	Plasma	Lui et al., 2014
Plasma	Bird	Gall et al., 2013
Rat	Plasma	Wu et al., 2015 ; Diwan et al., 2014 ; Choi et al., 2013 ; Tinwell et al., 2013
Rat	Serum	Jugdaohsingh et al., 2015 ; Strakovsky et al., 2014 ; Betz et al., 2013
Rat	Tissue Lysate	Zhang et al., 2014

For Technical Support:

techserv-usa@enzolifesciences.com

techserv-eu@enzolifesciences.com

Oxytocin ELISA Kit (ADI-900-153A)

SPECIES	SAMPLE ORIGIN	REFERENCES
Cat	Plasma	Bienboire-Frosini et al., 2017
Cow	Plasma	Bienboire-Frosini et al., 2017
Dog	Plasma	Bienboire-Frosini et al.;
Dog	Serum	Hollinshead et al., 2010
Goat	Plasma	Bienboire-Frosini et al., 2017
Horse	Plasma	Bienboire-Frosini et al., 2017
Human	Plasma	Sano et al., 2016 ; Baskaran et al., 2016 ; Husarova et al., 2016 ; Lin et al., 2015 ; Muin et al., 2015 ; Eidelman-Rothman et al., 2015 ; Emeny et al., 2015 ; Holt-Lunstad et al., 2008
Human	Saliva	Ormsby et al., 2016 ; Tsuji et al., 2015 ; Lebowitz et al., 2015 ; Jaeggi et al., 2015 ; Rubin et al., 2013 ; Dai et al., 2012 ; Feldman et al., 2010 ; Ray et al., 2009 ; Holt-Lunstad et al., 2008
Human	Urine	Cohen-Bendahan et al., 2015 ; Gregary et al., 2015 ; Salvante et al., 2012 ; Bick and Dozier, 2010
Human	CSF	Clark et al., 2013
Monkey	Urine	Finkenwirth et al., 2016 ; Finkenwirth et al., 2015 ; Moscovice and Zeigler, 2012 ; Snowdon et al., 2010
Monkey	Plasma	Freeman et al., 2016
Monkey	CSF	Freeman et al., 2016
Monkey	Serum	Michopoulos et al., 2011
Mouse	Plasma	Scott et al., 2015 ; Lopatina et al., 2011 ; Roizen et al., 2007
Pig	Plasma	Bienboire-Frosini et al., 2017 ; Norrby et al., 2011
Rabbit	Plasma	Szeto et al., 2013
Rabbit	Urine	Szeto et al., 2013
Rat	Plasma	Pereira, et al., 2015 ; Carini and Nephew, 2013 ; Vega et al., 2010 ; Martinez-Lorenzana et al., 2008
Rat	Tissue Lysate	Zhang et al., 2015
Rat	CSF	Martinez-Lorenzana et al., 2008
Seal	Plasma	Robinson et al., 2015
Sheep	Plasma	Bienboire-Frosini et al., 2017
Voles	Plasma	Yee et al., 2016 ; Sun et al., 2014

Progesterone ELISA Kit (ADI-900-011)

SPECIES	SAMPLE ORIGIN	REFERENCES
Cow	Serum	Mutinati et al., 2013
Cow	Plasma	Pothmann et al., 2015
Dog	Serum	Gabriel et al., 2016
Dolphin	Blubber	Kellar et al., 2017
Fish	Plasma	O'Connell et al., 2013 ; O'Connell and Hofmann, 2012
Horse	Plasma	Kuhl et al., 2016 ; Gajos et al., 2015
Horse	Serum	Siemieniuch et al., 2016
Horse	Saliva	Aurich et al., 2015
Human	Serum	Pietropolli et al., 2015
Human	Culture Supernatant	Ophir et al., 2014 ; Yung et al., 2014
Mouse	Plasma	Tanaka et al., 2016
Mouse	Culture Supernatant	Kim et al., 2016
Mouse	Serum	Facciponte et al., 2014
Octopus	Feces	Larson and Anderson, 2010
Pig	Serum	Dalmau et al., 2015
Rat	Plasma	Pyter et al., 2013 ; Tinwell et al., 2013
	Water	Kidd et al., 2013 ; Kidd et al., 2010

Testosterone ELISA Kit (ADI-900-065)

SPECIES	SAMPLE ORIGIN	REFERENCES
Bird	Plasma	Watts et al., 2016 ; Deviche et al., 2016 ; Davies et al., 2015 ; Lynn et al., 2015 ; Deviche et al., 2014 ; Watts and Hahn, 2012 ; DeVries et al., 2011
Fish	Serum	Singh and Srivastava, 2015
Fish	Plasma	Sessa et al., 2013 ; O'Connell et al., 2012 ; Greenwood et al., 2008
Gazelle	Feces	Ezenwa et al., 2012
Hamster	Serum	Greives et al., 2008 ; Demas et al., 2004
Hamster	Plasma	Wommack et al., 2004
Lizard	Plasma	Seddon and Hews, 2016
Mouse	Serum	Latsko et al., 2016 ; Thompson et al., 2015 ; de Peyster et al., 2008 ; Chang et al., 2004
Mouse	Culture Supernatant	Kim et al., 2016 ; Park et al., 2013
Mouse	Plasma	Volden et al., 2013 ; Manna et al., 2008
Mouse	Tissue Lysate	Ko et al., 2002
Naked mole rat	Serum	Peragine et al., 2016 ; Swift-Gallant et al., 2015
Octopus	Feces	Larson and Anderson, 2010
Rabbit	Plasma	Traish et al., 2003
Rat	Serum	Cyrenne and Brown, 2011
Rat	Plasma	Brouard et al., 2016 ; Diwan et al., 2014 ; Choi et al., 2013 ; Borrow et al., 2013
Rat	Rat hair	Arnon et al., 2016
Sea turtle	Plasma	Allen et al., 2015
	Water	Kidd et al., 2013 ; Kidd et al., 2010



For more information, go to:

www.enzolifesciences.com/wildlife-endocrinology