

Aldosterone ELISA Kit (ADI-900-173)

SPECIES	SAMPLE ORIGIN	REFERENCES
Dog	Plasma	Wang et al., 2015
Human	Plasma	Neal et al., 2016 ; Mears et al., 2016 ; Conway et al., 2014
Mouse	Urine	Zhang et al., 2015
Mouse	Plasma	Li et al., 2018 ; Hathaway et al., 2015 ; Qiao et al., 2014
Mouse	Tissue Lysate	Li et al., 2018
Mouse	Serum	Mittag et al., 2013
Rat	Serum	Nagao et al., 2019 ; Owagboriaye et al., 2018 ; Sakata et al., 2015 ; Singh et al., 2012
Rat	Plasma	Zhou et al., 2018 ; de Souza et al., 2018 ; Feldhaus et al., 2017 ; Hohl et al., 2016 ; More et al., 2016 ; Dhar et al., 2014
Rat	Urine	Ilatovskaya et al., 2019 ; Mansuri et al., 2017 ; Katatayama et al., 2013
Rat	Tissue Lysate	Kobayashi et al., 2017

Angiotensin II ELISA Kit (ADI-900-204)

SPECIES	SAMPLE ORIGIN	REFERENCES
Human	Tissue Lysate	Holappa et al., 2015
Human	Plasma	McDonald et al., 2018 ; Pooja et al., 2018
Mouse	Tissue Lysate	Li et al., 2018 ; Werner et al., 2016
Mouse	Plasma	Dougherty et al., 2019 ; Li et al., 2018 ; Chen et al., 2014 ; Tsukuda et al., 2012
Mouse	Serum	Mittag et al., 2013
Pig	Plasma	Jakob et al., 2012
Rabbit	Plasma	Ning et al., 2018
Rat	Tissue Lysate	Xu et al., 2016
Rat	Plasma	Fu et al., 2019 ; Zeng et al., 2019 ; Zhou et al., 2018 ; Hohl et al., 2016 ; Gao et al., 2016 ; Waghe et al., 2015 ; Park et al., 2015

Complete citation can be found on product's webpage

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

SPECIES	SAMPLE ORIGIN	REFERENCES
Dog	Saliva	MacLean et al., 2018
Human	Blood	Brondino et al., 2018
Human	Plasma	Oztan et al., 2017 ; Corbett et al., 2016 ; Carson et al., 2015 ; Idu et al., 2015
Human	Saliva	Weisman et al., 2013
Human	Serum	Bonavia et al., 2019 ; Rubin et al., 2017 ; Mears et al., 2016 ; Rubin et al., 2013 ; Dai et al., 2012
Human	Urine	Cohen-Bendahan et al., 2015 ; Gray et al., 2007
Monkey	Plasma	Larke et al., 2016
Mouse	Urine	Wang et al., 2019 ; Thelen et al., 2015 ; Zhang et al., 2015 ; Hopp et al., 2015 ; Zhang et al., 2012
Mouse	Serum	Kaneko et al., 2011
Mouse	Plasma	Thelen et al., 2015 ; Sauzeau et al., 2007
Rat	Plasma	Sims et al., 2017 ; More et al., 2016 ; 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	Greenwood et al., 2018 ; Sims et al., 2017 ; Zhang et al., 2015 ; Cuadra et al., 2014 ; Greenwood et al., 2014
Rat	Serum	Ren et al., 2015
Rat	Culture Supernatant	Knott et al., 2008
Vole	Plasma	Sun et al., 2014

Corticosterone ELISA Kit (ADI-900-097)

SPECIES	SAMPLE ORIGIN	REFERENCES
Alligator	Plasma	Finger et al., 2019 ; Hamilton et al., 2018 ; Finger et al., 2018
Bird	Plasma	Sarpong et al., 2019 ; Guindre-Parker and Rubenstein, 2019 ; Cooper et al., 2019 ; Jimeno et al., 2018 ; Newberry et al., 2018 ; Levin et al., 2018 ; Eikenaar et al., 2018 ; Gao and Deviche, 2018 ; Jimeno et al., 2018 ; Robert et al., 2018 ; Davies et al., 2018 ; Aharon-Rotman et al., 2017 ; Squire et al., 2017 ; Davies et al., 2017 ; Sorenson, 2016 ; Hennin et al., 2016 ; Ernst et al., 2016 ; Deviche et al., 2016 ; Madliger et al., 2015 ; Eikenaar et al., 2015 ; Huth and Archer, 2015 ; Lynn et al., 2015 ; Buttemer et al., 2015 ; Steenweg et al., 2015 ; Archer, 2015 ; Deviche et al., 2014 ; Liu and Swanson, 2014 ; Mark and Rubenstein, 2013 ; Love et al., 2005
Bird	Blood	Crino et al., 2018 ; Crino et al., 2017 ; Marasco et al., 2015
Bird	Tissue Lysate	Harris et al., 2016
Bullfrog Tadpole	Blood	Wilkens et al., 2019
Chameleon	Tissue Lysate	Matas et al., 2016
Chicken	Blood	Hedlund et al., 2019
Chicken	Plasma	Gualtieri et al., 2019 ; Nelson et al., 2018 ; Nazar et al., 2017 ; Anjum et al., 2016 ; Ericsson and Jensen, 2016 ; Ericsson et al., 2016 ; Furukawa et al., 2016 ; Tachibana et al., 2007
Chicken	Serum	Gurung et al., 2018 ; Gurung et al., 2018 ; Kolodziejski et al., 2018 ; Olanrewaju et al., 2016
Fish	Plasma	O'Connell et al., 2013
Human	Plasma	Xia et al., 2018
Human	Culture Supernatant	Slominski et al., 2005
Lizard	Plasma	Seddon and Hews, 2016
Lynx	Feces	Fanson et al., 2012
Mouse	Serum	Aubry et al., 2019 ; Baazim et al., 2019 ; Ince et al., 2019 ; Lundberg et al., 2019 ; Luijten et al., 2019 ; Dos Santos et al., 2019 ; Shah et al., 2019 ; Moreno et al., 2018 ; Woo et al., 2018 ; Seo et al., 2018 ; Collison et al., 2018 ; Quinn et al., 2018 ; Kinouchi et al., 2018 ; Caratti et al., 2018 ; Choi et al., 2017 ; Abulseoud et al., 2017 ; Aghajani et al., 2017 ; Lee et al., 2016 ; Socala et al., 2016 ; Abelson et al., 2016 ; Grimley et al., 2016 ; Latsko et al., 2016 ; Li et al., 2016 ; Cawthorn et al., 2016 ; Penney et al., 2016 ; Saito et al., 2016 ; Shushimita et al., 2016 ; Wong et al., 2015 ; Ghilan et al., 2015 ; Desbonnet et al., 2015 ; Yokota et al., 2015 ; Mishima et al., 2015 ; Zhang et al., 2015 ; Wang et al., 2015 ; Burkus et al., 2015 ; Liu et al., 2015 ; Sial et al., 2015 ; Yang et al., 2014 ; Aghajani et al., 2013 ; Siemen et al., 2011 ; Fernández et al., 2008 ; Thompson et al., 2008

Mouse	Plasma	Abdurakhmanova et al., 2019 ; Zhao et al., 2019 ; Li et al., 2019 ; Gorny et al., 2019 ; Sullivan et al., 2019 ; Niraula et al., 2018 ; Pyter et al., 2018 ; Timper et al., 2018 ; Station et al., 2018 ; Corder et al., 2018 ; Lee et al., 2018 ; Cabellero-Hernandez et al., 2017 ; Mayerhofer et al., 2017 ; Soualeh et al., 2017 ; Picard et al., 2016 ; Arase et al., 2016 ; Ramirez et al., 2016 ; Cabral et al., 2016 ; Castellano et al., 2016 ; Chavan et al., 2016 ; Chang 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 ; Rodriguez et al., 2015 ; Lara-Padilla et al., 2015 ; Uzturk et al., 2015 ; Yamada et al., 2015 ; Lin and Sibille, 2015 ; Liu et al., 2015 ; Vega-Rivera et al., 2015 ; Liu et al., 2015 ; Wei et al., 2014 ; Brooks et al., 2014 ; Volden et al., 2013 ; Philbert et al., 2012 ; Zhai et al., 2009
Mouse	Tissue Lysate	Frank et al., 2019 ; Mahapatra et al., 2005
Mouse	Culture Supernatant	Bae et al., 2016
Mouse	Amniotic Fluid	Xiang et al., 2016
Mouse	Urine	Hutchinson et al., 2012 ; Fitchett et al., 2005
Mouse	Feces	Siemen et al., 2011
Mouse	Blood	Mancuso et al., 2006
Penguin	Plasma	Stier et al., 2019
Pig	Culture Supernatant	Hu et al., 2017
Rat	Plasma	Strzelewicz et al., 2019 ; Bekhbat et al., 2019 ; Huzard et al., 2019 ; Mandal et al., 2019 ; Ali et al., 2019 ; Hayes et al., 2019 ; Hasan Mahmood et al., 2019 ; Ibrahim et al., 2019 ; Pitcher et al., 2019 ; Conoscenti et al., 2019 ; Satta et al., 2018 ; Huzard et al., 2018 ; Bekhbat et al., 2018 ; Kentner et al., 2018 ; Vore et al., 2017 ; Feldhaus et al., 2017 ; Katahira et al., 2017 ; Wang et al., 2017 ; Bülbül et al., 2017 ; Contreras et al., 2017 ; Torcaso et al., 2016 ; Bray et al., 2016 ; Bekhbat et al., 2016 ; Baracz et al., 2016 ; Karanges et al., 2016 ; Lee et al., 2016 ; Jones et al., 2016 ; Cason et al., 2016 ; Page et al., 2016 ; Gano et al., 2016 ; Patarroyo et al., 2016 ; Leo et al., 2015 ; Er et al., 2015 ; Ruvalcaba-Delgadillo 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 ; Wang et al., 2015 ; Khalil, 2015 ; Lajud et al., 2012 ; Charles et al., 2012 ; Aydin et al., 2011 ; Ray 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	Serum	Navarro-Meza et al., 2019 ; Chakraborty and Chattarji, 2019 ; Pan et al., 2019 ; Zhao et al., 2019 ; Owagboriaye et al., 2018 ; Fonken et al., 2018 ; Sun et al., 2018 ; Zhang et al., 2017 ; Kinn Rød et al., 2017 ; Bonini et al., 2016 ; Lin et al., 2016 ; Choi et al., 2016 ; Novaes et al., 2016 ; Williamson et al., 2016 ; Mou et al., 2016 ; Koe et al., 2016 ; Gano et al., 2016 ; Jacinto et al., 2016 ; Licznerski et al., 2015 ; Jing et al., 2015 ; Schneider et al., 2015 ; Nowland et al., 2011 ; Johnston et al., 2007
Rat	Feces	Sullivan et al., 2019 ; Green et al., 2018 ; Vega-Torres et al., 2018 ; Turner et al., 2012
Rat	Culture Supernatant	Sobesky et al., 2016 ; Charles et al., 2015
Rat	Hair	Arnon et al., 2016
Rat	Urine	Wang et al., 2011
Toad	Plasma	Gardner et al., 2018
Turtle	Plasma	West et al., 2018
Vole	Plasma	Soto et al., 2019 ; Jarcho et al., 2019 ; McNeal et al., 2018 ; Yee et al., 2016 ; Scotti et al., 2015

Cortisol ELISA Kit (ADI-900-071)

SPECIES	SAMPLE ORIGIN	REFERENCES
Cow	Plasma	Kelly et al., 2017 ; Olson et al., 2016 ; Earley et al., 2015
Fish	Serum	Khan et al., 2019
Fish	Culture Supernatant	Raja et al., 2019
Fish	Plasma	Wei et al., 2018 ; Kim and Kang, 2016 ; Kim and Kang, 2016 ; Kim and Kang, 2016 ; Delfosse et al., 2016 ; Mommer and Bell, 2013 ; Liss et al., 2013 ; Tapia et al., 2012 ; Greenwood et al., 2008
Fish	Tissue Lysate	King et al., 2016
Gazelle	Feces	Ezenwa et al., 2012
Hamster	Serum	Carlton and Demas, 2017 ; Demas et al., 2004 ; Wommack and Delville, 2003
Hamster	Plasma	Wommack et al., 2004
Horse	Plasma	Lemasson et al., 2015
Human	Plasma	Skornyakov et al., 2019 ; Watt et al., 2016
Human	Serum	Carrico et al., 2018
Human	Breast Milk	O'Rourke et al., 2018
Human	Saliva	Walton et al., 2016 ; Allen et al., 2016 ; Starnino et al., 2016 ; Nelson et al., 2008

Human	Culture Supernatant	Escajadillo et al., 2016 ; Lucki et al., 2012
Human	Urine	Fukuda et al., 2008
Lion	Feces	Creel et al., 2013
Monkey	Feces	Cinque et al., 2017
Mouse	Culture Supernatant	Tomabechi et al., 2018
Mouse	Serum	Han et al., 2018 ; Cawthorn et al., 2016 ; Gabaglia et al., 2007
Mouse	Plasma	Verma et al., 2016
Mouse	Tissue Lysate	Oh et al., 2015
Mouse	Feces	Running, 2015
Pig	Amniotic Fluid	Trus et al., 2018
Pig	Plasma	Trus et al., 2018 ; Turpin et al., 2016 ; Turpin et al., 2016 ; Rault et al., 2015
Pig	Saliva	Zupan et al., 2017
Rat	Amniotic Fluid	Jang et al., 2018
Rat	Plasma	Mehus et al., 2017
Sheep	Saliva	Häger et al., 2017
Whale	Culture Supernatant	Trumble et al., 2018
Whale	Breath Condensate	Thompson, et al., 2014
-	Water	Henderson et al., 2017

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	Plasma	Kozakiewicz et al., 2019 ; Duffy et al., 2015 ; Kumar et al., 2015 ; Li et al., 2014 ; Jeanes et al., 2004
Human	Tissue Lysate	Johnston et al., 2016
Human	Serum	Mukhtar et al., 2016 ; Puthuachary and Nathan, 2008
Human	Culture Supernatant	Patel and Bhat, 2004
Human	Urine	Huang et al., 2002
Mouse	Plasma	Hoffman-Goetz and Quadrilatero, 2003
Rat	Plasma	Goharinia et al., 2017
Rat	Tissue Lysate	Fang and Lin, 2008

DHEA ELISA Kit (ADI-900-093)

SPECIES	SAMPLE ORIGIN	REFERENCES
Cow	Serum	O'Driscoll et al., 2017
Cow	Plasma	O'Driscoll et al., 2015
Human	Tissue Lysate	Hong et al., 2019 ; Shin et al., 2018
Human	Serum	Hong et al., 2019 ; Chen et al., 2018 ; Shin et al., 2018 ; Huang et al., 2017 ; Sato et al., 2016
Human	Culture Supernatant	Zhu et al., 2019 ; Pyun et al., 2018 ; Chuang et al., 2017 ; Escajadillo et al., 2016 ; Lucki et al., 2012
Human	Plasma	Chuang et al., 2017
Rat	Feces	Sullivan et al., 2019
Rat	Culture Supernatant	Yang et al., 2018
Rat	Plasma	Horii et al., 2016

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

SPECIES	SAMPLE ORIGIN	REFERENCES
Fish	Culture Supernatant	Truong et al., 2016
Fish	Plasma	Sessa et al., 2013
Human	Serum	Mamun et al., 2018 ; Toonen et al., 2017
Human	Culture Supernatant	Xu et al., 2014 ; Bukovsky, 2009
Human	Cell Lysate	Fang et al., 2013
Mouse	Plasma	Zhang et al., 2019 ; Ear et al., 2019
Mouse	Culture Supernatant	Kim et al., 2018 ; Wang et al., 2014
Mouse	Serum	Chacraborty et al., 2016 ; Facciponte et al., 2014
Mouse	Tissue Lysate	Maurya et al., 2014
Octopus	Feces	Larson and Anderson, 2010
Rat	Plasma	Ear et al., 2019
Rat	Serum	Oludare et al., 2016
Rat	Culture Supernatant	Sittajody et al., 2013 ; Wang et al., 2013
Sea otter	Feces	Larson et al., 2013
-	Water	Heub et al., 2015 ; Kidd et al., 2013 ; Kidd et al., 2010 ; Hanselman et al., 2004

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

SPECIES	SAMPLE ORIGIN	REFERENCES
Bird	Blood	Crino et al., 2018
Bird	Plasma	Wilcoxon et al., 2015 ; Gall et al., 2013
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	Liu et al., 2014
Pig	Follicular Fluid	Pawlak et al., 2018
Rat	Plasma	Sittajody et al., 2017 ; Saiku and Vyas, 2017 ; Contreras and García, 2017 ; Tomaz et al., 2016 ; Wu et al., 2015 ; Diwan et al., 2014 ; Choi et al., 2013 ; Tinwell et al., 2013
Rat	Serum	Vieyra et al., 2019 ; Lee et al., 2016 ; Jugdaohsingh et al., 2015 ; Strakovsky et al., 2014 ; Betz et al., 2013 ; Tseng et al., 2012
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
Cow	Saliva	Geburt et al., 2015
Dog	Saliva	MacLean et al., 2018
Dog	Plasma	Petersson et al., 2017
Dog	Serum	Hollinshead et al., 2010
Horse	Plasma	Lansade et al., 2018 ; Malinkowski et al., 2018 ; Ginther and Beg, 2011
Human	Plasma	Hayashi et al., 2019 ; Kampka et al., 2019 ; Brondino et al., 2018 ; Oztan et al., 2017 ; Sano et al., 2016 ; Corbett et al., 2016 ; Kosaka et al., 2016 ; Husarova et al., 2016 ; Lin et al., 2015 ; Muin et al., 2015 ; Eidelman-Rothman et al., 2015 ; Emeny et al., 2015 ; Szeto et al., 2011 ; Holt-Lunstad et al., 2008 ; Morhenn et al., 2008
Human	Saliva	Takahata et al., 2018 ; Behr et al., 2017 ; Ormsby et al., 2016 ; Tsuji et al., 2015 ; Lebowitz et al., 2015 ; Jaeggi et al., 2015 ; Feldman et al., 2010 ; Ray et al., 2009 ; Holt-Lunstad et al., 2008
Human	Serum	Miwa et al., 2018 ; Behr et al., 2017 ; Rubin et al., 2017 ; Sasaki et al., 2016 ; Lin et al., 2013 ; Rubin et al., 2013 ; Dai et al., 2012 ; Morhenn et al., 2008
Human	Urine	Eisenberg et al., 2018 ; Algoe et al., 2017 ; Cohen-Bendahan et al., 2015 ; Gregory 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 Ziegler, 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	Ear et al., 2019 ; Damoah et al., 2019 ; Denizot et al., 2016 ; Larke et al., 2016 ; Scott et al., 2015 ; Lopatina et al., 2011 ; Roizen et al., 2007
Mouse	CSF	Damoah et al., 2019
Pig	Plasma	Rius et al., 2018 ; Norrby et al., 2011
Rabbit	Plasma	Szeto et al., 2013
Rat	Plasma	Holubová et al., 2019 ; Ear et al., 2019 ; Baracz et al., 2016 ; Pereira et al., 2015 ; Carini and Nephew, 2013 ; Vega et al., 2010 ; Martínez-Lorenzana et al., 2008
Rat	Serum	Liberto et al., 2018 ; Inoue et al., 2016
Rat	Tissue Lysate	Meidahl et al., 2018 ; Zhang et al., 2015 ; Martínez-Lorenzana et al., 2008
Rat	Culture Supernatant	Maejima et al., 2017
Rat	CSF	Couto-Pereira et al., 2016 ; Martínez-Lorenzana et al., 2008
Rat	Microdialysates	Bülbul et al., 2016
Seal	Plasma	Robinson et al., 2015
Vole	Plasma	Yee et al., 2016 ; Sun et al., 2014

For Technical Support:
 techserv-usa@enzolifesciences.com
 techserv-eu@enzolifesciences.com

Progesterone ELISA Kit (ADI-900-011)

SPECIES	SAMPLE ORIGIN	REFERENCES
Bear	Hair	Cattet et al., 2018
Cat	Culture Supernatant	Jursza-Piotrowska et al., 2016 ; Siemieniuch et al., 2014
Cow	Culture Supernatant	Alemu et al., 2018
Cow	Plasma	Wagener et al., 2017 ; Andreas et al., 2016 ; Pothmann et al., 2015
Cow	Follicular Fluid	Rizzo et al., 2016
Cow	Serum	Ahmed et al., 2016 ; Mutinati et al., 2013
Dog	Serum	Gabriel et al., 2016
Dolphin	Blubber	Trego et al., 2019
Fish	Plasma	O'Connell et al., 2013 ; O'Connell and Hofmann, 2012 ; Kidd et al., 2010
Horse	Plasma	Aurich et al., 2019 ; Melchert et al., 2019 ; Kuhl et al., 2016 ; Gajos et al., 2015 ; Köhne et al., 2014
Horse	Endometrial Fluid	Beyer et al., 2018
Horse	Serum	Siemieniuch et al., 2017 ; Köhne et al., 2016 ; Siemieniuch et al., 2016
Horse	Follicular Fluid	Scarlet et al., 2015
Horse	Saliva	Aurich et al., 2015
Human	Follicular Fluid	O'Brien et al., 2019
Human	Culture Supernatant	Sagers et al., 2018 ; Ophir et al., 2014 ; Yung et al., 2014
Human	Serum	Pietropoli et al., 2015
Mouse	Plasma	Ishii et al., 2019 ; Tanaka et al., 2016
Mouse	Serum	Nakano et al., 2018 ; Ziegler-Waldkirch et al., 2018 ; Filipovich et al., 2016 ; Facciponte et al., 2014
Mouse	Culture Supernatant	Mittelman-Smith et al., 2018 ; Kim et al., 2016
Octopus	Feces	Larson and Anderson, 2010
Pig	Follicular Fluid	Pawlak et al., 2018
Pig	Serum	Dalmau et al., 2015
Rat	Serum	Vieyra et al., 2019 ; Oludare et al., 2016
Rat	Culture Supernatant	Elmorsy et al., 2017 ; Qiu et al., 2017 ; Sittajody et al., 2013 ; Wang et al., 2013
Rat	Plasma	Sittajody et al., 2017 ; Saiku and Vyas, 2017 ; Contreras and Gutiérrez-García, 2017 ; Pyter et al., 2013 ; Tinwell et al., 2013
Whale	Blubber	Pallin et al., 2018 ; Kellar et al., 2013
Whale	Serum	Kellar et al., 2013
Whale	Urine	Kellar et al., 2013
-	Water	Kidd et al., 2013 ; Kidd et al., 2010

For Technical Support:
techserv-usa@enzolifesciences.com
techserv-eu@enzolifesciences.com

Testosterone ELISA Kit (ADI-900-065)

SPECIES	SAMPLE ORIGIN	REFERENCES
Bat	Serum	Souza et al., 2018
Bear	Hair	Cattet et al., 2018
Bird	Plasma	Graham et al., 2019 ; Levin et al., 2018 ; Robart et al., 2018 ; Davies et al., 2018 ; Needham et al., 2017 ; Li et al., 2017 ; Whittaker et al., 2017 ; Watts et al., 2016 ; Alward et al., 2016 ; Ernst et al., 2016 ; Deviche et al., 2016 ; Bauer et al., 2016 ; Sorensen et al., 2016 ; Davies et al., 2015 ; Wilcoxon et al., 2015 ; Lynn et al., 2015 ; Davies et al., 2015 ; Deviche et al., 2014 ; Rosvall, 2013 ; Watts and Hahn, 2012 ; DeVries et al., 2011
Bird	Serum	Alward et al., 2016
Bird	Dialysate	Ikeda et al., 2014
Dog	Culture Supernatant	Lea et al., 2016
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
Gerbil	Serum	Zanatelli et al., 2014
Hamster	Serum	Long et al., 2018 ; Greives et al., 2008 ; Demas et al., 2004
Hamster	Plasma	Wommack et al., 2004
Horse	Plasma	Lemasson et al., 2015
Human	Serum	Bawor et al., 2014
Lizard	Plasma	LaDage et al., 2017 ; Seddon and Hews, 2016
Mouse	Serum	Templin et al., 2019 ; Weger et al., 2019 ; Hu et al., 2018 ; Mamun et al., 2018 ; Lee et al., 2018 ; Zhang et al., 2017 ; Toonen et al., 2017 ; Latsko et al., 2016 ; Thompson et al., 2015 ; de Peyster et al., 2008 ; Chang et al., 2004
Mouse	Plasma	Shawki et al., 2018 ; Fu et al., 2018 ; Lopes and König, 2016 ; Homma et al., 2016 ; Volden et al., 2013 ; Manna et al., 2008
Mouse	Culture Supernatant	Zhang et al., 2017 ; Kim et al., 2016 ; Park et al., 2013
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	Souza et al., 2019 ; Ribeiro et al., 2019 ; Migliaccio et al., 2019 ; Ribeiro et al., 2018 ; Moon et al., 2017 ; Zhu et al., 2017 ; Lynch 3rd et al., 2016 ; Ha et al., 2016 ; Choi et al., 2016 ; Lynch 3rd et al., 2016 ; Jin et al., 2013 ; Tseng et al., 2012 ; Cyrenne and Brown, 2011
Rat	Plasma	Asimes et al., 2019 ; Sittajody et al., 2017 ; Saiku and Vyas, 2017 ; Brouard et al., 2016 ; Yamaguchi et al., 2016 ; Thangapandian and Miltonprabu, 2015 ; Diwan et al., 2014 ; Choi et al., 2013 ; Borrow et al., 2013 ; Tinwell et al., 2013
Rat	Hair	Arnon et al., 2016
Rat	Tissue Lysate	Clarke and Pearl, 2014
Snail	Hemolymph	Ibrahim and Sayed, 2019
Snail	Tissue Lysate	Ibrahim and Ghoname, 2018 ; Omran and Salama, 2016
Turtle	Plasma	Jensen et al., 2018 ; Allen et al., 2015
Whale	Blubber	Cates et al., 2019
-	Water	Kidd et al., 2013 ; Kidd et al., 2010

For more information, go to:

www.enzolifesciences.com/wildlife-endocrinology