

## GRO-CINC-1 (rat) ELISA Kit (ADI-900-074)

For the quantitative determination of rat GRO/CINC-1 in culture supernatants, plasma, and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
BALF	Rat	G.T. Wagenaar, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24951776">https://www.ncbi.nlm.nih.gov/pubmed/24951776</a>
BALF	Rat	X. Chen, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26495902">https://www.ncbi.nlm.nih.gov/pubmed/26495902</a>
Intestinal tissue	Rat	V.R. Yadav, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27503458">https://www.ncbi.nlm.nih.gov/pubmed/27503458</a>
Liver tissue	Rat	V.R. Yadav, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27503458">https://www.ncbi.nlm.nih.gov/pubmed/27503458</a>
Lung tissue	Rat	K. Hashimoto, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26725713">https://www.ncbi.nlm.nih.gov/pubmed/26725713</a>
Plasma	Rat	V.R. Yadav, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/25645333">https://www.ncbi.nlm.nih.gov/pubmed/25645333</a>
Plasma	Rat	V.R. Yadav, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27503458">https://www.ncbi.nlm.nih.gov/pubmed/27503458</a>

## IFN- $\gamma$ (human) ELISA Kit (ADI-900-136)

For the quantitative determination of human IFN- $\gamma$  in culture supernatants, plasma, serum, and urine.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Culture supernatant	Human	L. Hu, <i>et al.</i> (2012)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22406996">https://www.ncbi.nlm.nih.gov/pubmed/22406996</a>
Culture supernatant	Human	R. Maitra, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28422714">https://www.ncbi.nlm.nih.gov/pubmed/28422714</a>
Saliva	Human	E.C. Siqueira, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26876491">https://www.ncbi.nlm.nih.gov/pubmed/26876491</a>

## IFN- $\gamma$ (mouse) ELISA Kit (ADI-900-137)

For the quantitative determination of mouse IFN- $\gamma$  in culture supernatants, plasma, serum, and urine.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Culture supernatant	Mouse	Y.S. Wu, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24723960">https://www.ncbi.nlm.nih.gov/pubmed/24723960</a>
Culture supernatant	Mouse	C.Y. Wang, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27520377">https://www.ncbi.nlm.nih.gov/pubmed/27520377</a>
Serum	Mouse	M.S. Kim, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24530874">https://www.ncbi.nlm.nih.gov/pubmed/24530874</a>
Serum	Mouse	Y.S. Wu, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24723960">https://www.ncbi.nlm.nih.gov/pubmed/24723960</a>
Serum	Mouse	Y. Aoyama, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/25886696">https://www.ncbi.nlm.nih.gov/pubmed/25886696</a>
Serum	Mouse	C.Y. Wang, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27520377">https://www.ncbi.nlm.nih.gov/pubmed/27520377</a>

Advance Your Research:  
[www.enzolifesciences.com](http://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

## IGF-1 (human) ELISA Kit (ADI-900-150)

For the quantitative determination of human IGF-1 in plasma and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Culture supernatant	Human	L.M. McGinley, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26744412">https://www.ncbi.nlm.nih.gov/pubmed/26744412</a>
Drug microparticles	Human	W. Lee, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27381562">https://www.ncbi.nlm.nih.gov/pubmed/27381562</a>
Plasma	Human	W.D. Arnold, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27907033">https://www.ncbi.nlm.nih.gov/pubmed/27907033</a>
Serum	Human	M. Spillane, <i>et al.</i> (2011)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22050827">https://www.ncbi.nlm.nih.gov/pubmed/22050827</a>
Serum	Human	W.R. Duff, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24281841">https://www.ncbi.nlm.nih.gov/pubmed/24281841</a>
Serum	Human	M. Spillane, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/25435783">https://www.ncbi.nlm.nih.gov/pubmed/25435783</a>
Serum	Human	T. Floros, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26732162">https://www.ncbi.nlm.nih.gov/pubmed/26732162</a>
Serum	Human	M. Spillane, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26957922">https://www.ncbi.nlm.nih.gov/pubmed/26957922</a>
Velvet antler	Deer	S.H. Tseng, <i>et al.</i> (2012)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22951396">https://www.ncbi.nlm.nih.gov/pubmed/22951396</a>

## IL-1 $\beta$ (human) ELISA Kit (ADI-900-130A)

For the quantitative determination of IL-1 $\beta$  in culture supernatant, plasma, serum, and urine of human origin.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Culture supernatant	Human	S.Y. Park, <i>et al.</i> (2010)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/20131233">https://www.ncbi.nlm.nih.gov/pubmed/20131233</a>
Culture supernatant	Human	S. Raha, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27477351">https://www.ncbi.nlm.nih.gov/pubmed/27477351</a>
Neocortical tissue	Human	L. Lorigados-Pedre, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29401729">https://www.ncbi.nlm.nih.gov/pubmed/29401729</a>
Saliva	Human	E.C. Siqueira, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26876491">https://www.ncbi.nlm.nih.gov/pubmed/26876491</a>
Serum	Human	L. Lorigados-Pedre, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29401729">https://www.ncbi.nlm.nih.gov/pubmed/29401729</a>

## IL-1 $\beta$ (rat) ELISA Kit (ADI-900-131A)

For the quantitative determination of IL-1 $\beta$  in culture supernatants and serum of rat origin.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Hippocampus	Rat	H. Fan, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23885269">https://www.ncbi.nlm.nih.gov/pubmed/23885269</a>

## IL-2 (human) ELISA Kit (ADI-900-118A)

For the quantitative determination of human IL-2 in culture supernatants, plasma, and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Serum	Human	G.U. Hong, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30195017">https://www.ncbi.nlm.nih.gov/pubmed/30195017</a>

## IL-2 (mouse) ELISA Kit (ADI-900-042)

For the quantitative determination of mouse IL-2 in culture supernatants and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Brain tissue	Mouse	G.U. Hong, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30195017">https://www.ncbi.nlm.nih.gov/pubmed/30195017</a>
Serum	Mouse	K. Yoshino, <i>et al.</i> (2008)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/18387131">https://www.ncbi.nlm.nih.gov/pubmed/18387131</a>
Serum	Mouse	G.U. Hong, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30195017">https://www.ncbi.nlm.nih.gov/pubmed/30195017</a>

## IL-4 (human) ELISA Kit (ADI-900-145A)

For the quantitative determination of human IL-4 in culture supernatants, plasma, and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Serum	Human	A. Milajerdi, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29641735">https://www.ncbi.nlm.nih.gov/pubmed/29641735</a>

## IL-4 (mouse) ELISA Kit (ADI-900-043)

For the quantitative determination of mouse IL-4 in culture supernatants and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Plasma	Mouse	Y. Zhao, <i>et al.</i> (2009)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/19156219">https://www.ncbi.nlm.nih.gov/pubmed/19156219</a>
Serum	Mouse	X. Ding, <i>et al.</i> (2019)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30587058">https://www.ncbi.nlm.nih.gov/pubmed/30587058</a>

## IL-6 (human) high sensitivity ELISA Kit (ENZ-KIT178)

For the highly sensitive quantification of human IL-6 in culture supernatants, plasma, serum, and urine.

SAMPLE TYPE	SPECIES	REFERENCES	LINK

## IL-6 (mouse) ELISA Kit (ADI-900-045)

For the quantitative determination of mouse IL-6 in culture supernatants and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Blood	Mouse	L. Cui, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27261584">https://www.ncbi.nlm.nih.gov/pubmed/27261584</a>
Brain tissue	Mouse	M. Taniguchi, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27892528">https://www.ncbi.nlm.nih.gov/pubmed/27892528</a>
Culture supernatant	Mouse	K. Lee, <i>et al.</i> (2012)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22578249">https://www.ncbi.nlm.nih.gov/pubmed/22578249</a>
Culture supernatant	Mouse	M.A. Hokenson, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23527226">https://www.ncbi.nlm.nih.gov/pubmed/23527226</a>
Culture supernatant	Mouse	Y.M. Shin, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23054007">https://www.ncbi.nlm.nih.gov/pubmed/23054007</a>
Culture supernatant	Mouse	P.S. Nayak, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26006045">https://www.ncbi.nlm.nih.gov/pubmed/26006045</a>
Culture supernatant	Mouse	J.Y. Seo, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26548345">https://www.ncbi.nlm.nih.gov/pubmed/26548345</a>
Culture supernatant	Mouse	Y.A. Bae, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27382358">https://www.ncbi.nlm.nih.gov/pubmed/27382358</a>
Culture supernatant	Mouse	J. Lee, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26709074">https://www.ncbi.nlm.nih.gov/pubmed/26709074</a>
Culture supernatant	Mouse	C.Y. Wang, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27520377">https://www.ncbi.nlm.nih.gov/pubmed/27520377</a>
Lung tissue	Mouse	A. Hsu, <i>et al.</i> (2007)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/17822444">https://www.ncbi.nlm.nih.gov/pubmed/17822444</a>
Plasma	Mouse	A.V. Aubry, <i>et al.</i> (2019)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30542090">https://www.ncbi.nlm.nih.gov/pubmed/30542090</a>
Serum	Mouse	S.Y. Park, <i>et al.</i> (2010)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/20131233">https://www.ncbi.nlm.nih.gov/pubmed/20131233</a>
Serum	Mouse	C.Y. Wang, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27520377">https://www.ncbi.nlm.nih.gov/pubmed/27520377</a>
Serum	Mouse	B. Shan, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28346409">https://www.ncbi.nlm.nih.gov/pubmed/28346409</a>

## IL-8 (human) ELISA Kit (ADI-900-156)

For the quantitative determination of human IL-8 in culture supernatants, plasma, and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Culture supernatant	Human	A.N. Smith, <i>et al.</i> (2012)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22105830">https://www.ncbi.nlm.nih.gov/pubmed/22105830</a>
Serum	Human	A.C. Nilsson, <i>et al.</i> (2008)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/18356328">https://www.ncbi.nlm.nih.gov/pubmed/18356328</a>
Serum	Human	C.D. Ene, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26504364">https://www.ncbi.nlm.nih.gov/pubmed/26504364</a>

## IL-10 (human) ELISA Kit (ADI-900-036)

For the quantitative determination of human IL-10 in culture supernatants, plasma, and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Saliva	Human	E.C. Siqueira, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26876491">https://www.ncbi.nlm.nih.gov/pubmed/26876491</a>

## IL-12p70 (human) ELISA Kit (ADI-900-202)

For the quantitative determination of human IL-12p70 in culture supernatants, plasma, and serum samples.

SAMPLE TYPE	SPECIES	REFERENCES	LINK

## IL-13 (human) ELISA Kit (ADI-900-208)

For the quantitative determination of human IL-13 in plasma, serum and cell supernatant samples.

SAMPLE TYPE	SPECIES	REFERENCES	LINK

### IL-17A (human) ELISA Kit (ADI-900-177)

For the quantitative determination of human IL-17A in culture supernatants, plasma, and serum samples.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Serum	Human	D. Golan, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23767916">https://www.ncbi.nlm.nih.gov/pubmed/23767916</a>

### IL-33 (human) ELISA Kit (ADI-900-201)

For the quantitative determination of human IL-33 in cell lysates, culture supernatants, serum, plasma and synovial fluid samples.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Cell lysate	Human	R. Kakkar, <i>et al.</i> (2012)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22215666">https://www.ncbi.nlm.nih.gov/pubmed/22215666</a>
Culture supernatant	Human	R. Kakkar, <i>et al.</i> (2012)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22215666">https://www.ncbi.nlm.nih.gov/pubmed/22215666</a>
Serum	Human	J. Yue, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27026387">https://www.ncbi.nlm.nih.gov/pubmed/27026387</a>

### Leptin (human) ELISA Kit (ADI-900-028A)

For the quantitative determination of human leptin in culture supernatants, plasma, and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Plasma	Human	Y.H. Lin, <i>et al.</i> (2011)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/21195702">https://www.ncbi.nlm.nih.gov/pubmed/21195702</a>
Plasma	Human	G. Derosa, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23467997">https://www.ncbi.nlm.nih.gov/pubmed/23467997</a>
Plasma	Human	G. Derosa, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/25994603">https://www.ncbi.nlm.nih.gov/pubmed/25994603</a>
Plasma	Human	M. Tellechea, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/25598199">https://www.ncbi.nlm.nih.gov/pubmed/25598199</a>
Plasma	Human	B. Almeida-Pititto, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30479668">https://www.ncbi.nlm.nih.gov/pubmed/30479668</a>
Serum	Human	E.K. Kim, <i>et al.</i> (2011)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/21745625">https://www.ncbi.nlm.nih.gov/pubmed/21745625</a>
Serum	Human	Y. Ma, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24167669">https://www.ncbi.nlm.nih.gov/pubmed/24167669</a>
Serum	Human	R.K. Joshi, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24976730">https://www.ncbi.nlm.nih.gov/pubmed/24976730</a>
Serum	Human	E. Chavarria-Avila, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26064921">https://www.ncbi.nlm.nih.gov/pubmed/26064921</a>
Serum	Human	R. Rehman, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29805405">https://www.ncbi.nlm.nih.gov/pubmed/29805405</a>

### Leptin (mouse) ELISA Kit (ADI-900-019A)

For the quantitative determination of mouse leptin in culture supernatants, plasma, and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Culture supernatant	Mouse	S.D. Vangipuram, <i>et al.</i> (2007)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/16703005">https://www.ncbi.nlm.nih.gov/pubmed/16703005</a>
Liver tissue	Mouse	S. Nepali, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29484428">https://www.ncbi.nlm.nih.gov/pubmed/29484428</a>
Lung tissue	Mouse	A. Hsu, <i>et al.</i> (2007)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/17822444">https://www.ncbi.nlm.nih.gov/pubmed/17822444</a>
Plasma	Mouse	J. Wang, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26361146">https://www.ncbi.nlm.nih.gov/pubmed/26361146</a>
Plasma	Mouse	P.R. Burghardt, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26926827">https://www.ncbi.nlm.nih.gov/pubmed/26926827</a>
Plasma	Mouse	S.C. Fordahl, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27267686">https://www.ncbi.nlm.nih.gov/pubmed/27267686</a>
Plasma	Mouse	A. Sassmann-Schweda, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27699262">https://www.ncbi.nlm.nih.gov/pubmed/27699262</a>
Plasma	Mouse	S.H. Hong, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29474103">https://www.ncbi.nlm.nih.gov/pubmed/29474103</a>
Plasma	Mouse	M. Woo, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30200239">https://www.ncbi.nlm.nih.gov/pubmed/30200239</a>
Serum	Mouse	G.K. Bhat, <i>et al.</i> (2003)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/12606383">https://www.ncbi.nlm.nih.gov/pubmed/12606383</a>
Serum	Mouse	M.L. Hamm, <i>et al.</i> (2004)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/14985253">https://www.ncbi.nlm.nih.gov/pubmed/14985253</a>
Serum	Mouse	L. Recinella, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24021480">https://www.ncbi.nlm.nih.gov/pubmed/24021480</a>
Serum	Mouse	H.N. Choi, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29018783">https://www.ncbi.nlm.nih.gov/pubmed/29018783</a>
Serum	Mouse	K.A. Romano, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28844887">https://www.ncbi.nlm.nih.gov/pubmed/28844887</a>
Serum	Mouse	B. Shan, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28346409">https://www.ncbi.nlm.nih.gov/pubmed/28346409</a>
Serum	Mouse	I.H. Kim, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29207025">https://www.ncbi.nlm.nih.gov/pubmed/29207025</a>

## Leptin (rat) ELISA Kit (ADI-900-015A)

For the quantitative determination of rat leptin in culture supernatants, plasma, and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Plasma	Rat	B. Merino, <i>et al.</i> (2008)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/18096657">https://www.ncbi.nlm.nih.gov/pubmed/18096657</a>
Plasma	Rat	B. Galvez-Prieto, <i>et al.</i> (2012)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22679436">https://www.ncbi.nlm.nih.gov/pubmed/22679436</a>
Plasma	Rat	M. Choi, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23548559">https://www.ncbi.nlm.nih.gov/pubmed/23548559</a>
Plasma	Rat	J.C. de Oliveira, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23151360">https://www.ncbi.nlm.nih.gov/pubmed/23151360</a>
Plasma	Rat	S.A. Messenger, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22364734">https://www.ncbi.nlm.nih.gov/pubmed/22364734</a>
Plasma	Rat	S.A. Messenger, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23201827">https://www.ncbi.nlm.nih.gov/pubmed/23201827</a>
Plasma	Rat	R. Barazzoni, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23512916">https://www.ncbi.nlm.nih.gov/pubmed/23512916</a>
Plasma	Rat	H.E. Yim, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23333086">https://www.ncbi.nlm.nih.gov/pubmed/23333086</a>
Plasma	Rat	T.H. Yang, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28911493">https://www.ncbi.nlm.nih.gov/pubmed/28911493</a>
Plasma	Rat	R. Mukherjee, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26116537">https://www.ncbi.nlm.nih.gov/pubmed/26116537</a>
Plasma	Rat	A. Malta, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27561682">https://www.ncbi.nlm.nih.gov/pubmed/27561682</a>
Plasma	Rat	J. Ciriello, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27222924">https://www.ncbi.nlm.nih.gov/pubmed/27222924</a>
Plasma	Rat	H.C. Liu, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28911640">https://www.ncbi.nlm.nih.gov/pubmed/28911640</a>
Plasma	Rat	M.M. Rahman, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28578257">https://www.ncbi.nlm.nih.gov/pubmed/28578257</a>
Serum	Rat	J.F. Ge, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23619134">https://www.ncbi.nlm.nih.gov/pubmed/23619134</a>
Serum	Rat	S. Li, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24489777">https://www.ncbi.nlm.nih.gov/pubmed/24489777</a>

## MCP-1 (rat) ELISA Kit (ADI-900-077)

For the quantitative determination of rat MCP-1 in culture supernatants and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Culture supernatant	Rat	K. Afrasiabi, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26124920">https://www.ncbi.nlm.nih.gov/pubmed/26124920</a>
Serum	Rat	S. Li, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24489777">https://www.ncbi.nlm.nih.gov/pubmed/24489777</a>
Serum	Rat	B. Shan, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28346409">https://www.ncbi.nlm.nih.gov/pubmed/28346409</a>

## Nampt (Visfatin/PBEF) (human) intracellular ELISA Kit (AG-45A-0006EK)

For the quantitative determination of human Nampt (visfatin/PBEF) in cell lysates.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Cell lysate	Human	C. Bian, <i>et al.</i> (2019)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30741937">https://www.ncbi.nlm.nih.gov/pubmed/30741937</a>
Culture supernatant	Human	T. Yonezawa, <i>et al.</i> (2006)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/17123517">https://www.ncbi.nlm.nih.gov/pubmed/17123517</a>
Culture supernatant	Human	A. Körner, <i>et al.</i> (2007)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/17878256">https://www.ncbi.nlm.nih.gov/pubmed/17878256</a>
Plasma	Human	E. Parvaresh Rizi, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26308293">https://www.ncbi.nlm.nih.gov/pubmed/26308293</a>
Plasma	Human	R.M. de Guia, <i>et al.</i> (2019)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/31207144">https://www.ncbi.nlm.nih.gov/pubmed/31207144</a>
Serum	Human	A. Körner, <i>et al.</i> (2007)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/17878256">https://www.ncbi.nlm.nih.gov/pubmed/17878256</a>
Serum	Human	M. Fasshauer, <i>et al.</i> (2008)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/18034779">https://www.ncbi.nlm.nih.gov/pubmed/18034779</a>
Serum	Human	R. Retnakaran, <i>et al.</i> (2008)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/18410550">https://www.ncbi.nlm.nih.gov/pubmed/18410550</a>
Serum	Human	C. Benedict, <i>et al.</i> (2012)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22090280">https://www.ncbi.nlm.nih.gov/pubmed/22090280</a>
Serum	Human	M. Nourbakhsh, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/25723377">https://www.ncbi.nlm.nih.gov/pubmed/25723377</a>



## Nampt (Visfatin/PBEF) (mouse/rat) dual ELISA Kit (AG-45A-0007EK)

For the quantitative determination of Nampt (Visfatin/PBEF) in mouse or rat serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Cell lysate	Mouse	J.S. Kim, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/25566462">https://www.ncbi.nlm.nih.gov/pubmed/25566462</a>
Cell lysate	Mouse	D.F. Pisani, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27386154">https://www.ncbi.nlm.nih.gov/pubmed/27386154</a>

## TGF- $\beta$ 1 ELISA Kit (ADI-900-155)

For the quantitative determination of TGF- $\beta$ 1 in culture supernatants, plasma, and serum of human, mouse, rat, and cow origin.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Aortic homogenate	Human	M.E. Soto, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27163200">https://www.ncbi.nlm.nih.gov/pubmed/27163200</a>
Aortic homogenate	Human	M.E. Soto, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29483877">https://www.ncbi.nlm.nih.gov/pubmed/29483877</a>
BALF	Human	Y. Liu, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28498479">https://www.ncbi.nlm.nih.gov/pubmed/28498479</a>
Brain tissue	Mouse	G.U. Hong, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30195017">https://www.ncbi.nlm.nih.gov/pubmed/30195017</a>
Cell lysate	Human	G. Longo, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27031101">https://www.ncbi.nlm.nih.gov/pubmed/27031101</a>
Culture supernatant	Human	L.C. Moraes, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26249205">https://www.ncbi.nlm.nih.gov/pubmed/26249205</a>
Culture supernatant	Human	Y. Liu, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28498479">https://www.ncbi.nlm.nih.gov/pubmed/28498479</a>
Culture supernatant	Human	S. Rosini, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29844053">https://www.ncbi.nlm.nih.gov/pubmed/29844053</a>
Culture supernatant	Mouse	M.G. Morales, <i>et al.</i> (2012)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22964022">https://www.ncbi.nlm.nih.gov/pubmed/22964022</a>
Culture supernatant	Mouse	S.H. Choi, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/25935484">https://www.ncbi.nlm.nih.gov/pubmed/25935484</a>
Culture supernatant	Rat	N. Frara, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28399671">https://www.ncbi.nlm.nih.gov/pubmed/28399671</a>
Liposomes	Human	C.B.M. Platania, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28973964">https://www.ncbi.nlm.nih.gov/pubmed/28973964</a>
Serum	Human	Y. Liu, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28498479">https://www.ncbi.nlm.nih.gov/pubmed/28498479</a>
Serum	Human	G.U. Hong, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30195017">https://www.ncbi.nlm.nih.gov/pubmed/30195017</a>
Serum	Mouse	G.U. Hong, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30195017">https://www.ncbi.nlm.nih.gov/pubmed/30195017</a>
Vitreous humor	Rabbit	C.B.M. Platania, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28973964">https://www.ncbi.nlm.nih.gov/pubmed/28973964</a>

## TNF- $\alpha$ (human) ELISA Kit (ADI-900-099)

For the quantitative determination of human TNF- $\alpha$  in culture supernatants, plasma, and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Culture supernatant	Human	S.Y. Park, <i>et al.</i> (2010)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/20131233">https://www.ncbi.nlm.nih.gov/pubmed/20131233</a>
Culture supernatant	Human	S. Raha, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27477351">https://www.ncbi.nlm.nih.gov/pubmed/27477351</a>
Culture supernatant	Human	C.B.M. Platania, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30013474">https://www.ncbi.nlm.nih.gov/pubmed/30013474</a>
Plasma	Human	K.A. Kim, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27812515">https://www.ncbi.nlm.nih.gov/pubmed/27812515</a>
Plasma	Human	F. Falasca, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28830393">https://www.ncbi.nlm.nih.gov/pubmed/28830393</a>
Serum	Human	S. Buscemi, <i>et al.</i> (2012)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/22492368">https://www.ncbi.nlm.nih.gov/pubmed/22492368</a>
Serum	Human	A. Rangel-Lopez, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23408844">https://www.ncbi.nlm.nih.gov/pubmed/23408844</a>
Serum	Human	R.K. Joshi, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24976730">https://www.ncbi.nlm.nih.gov/pubmed/24976730</a>
Serum	Human	H.N. Choi, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29018783">https://www.ncbi.nlm.nih.gov/pubmed/29018783</a>
Serum	Human	G.U. Hong, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30195017">https://www.ncbi.nlm.nih.gov/pubmed/30195017</a>
Serum	Human	A. Milajerdi, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29641735">https://www.ncbi.nlm.nih.gov/pubmed/29641735</a>

## TNF- $\alpha$ (mouse) ELISA kit (ADI-900-047)

For the quantitative determination of human TNF- $\alpha$  in culture supernatants, plasma, and serum.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Blood	Mouse	L. Cui, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27261584">https://www.ncbi.nlm.nih.gov/pubmed/27261584</a>
Brain tissue	Mouse	G.U. Hong, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30195017">https://www.ncbi.nlm.nih.gov/pubmed/30195017</a>
Culture supernatant	Mouse	Y.M. Shin, <i>et al.</i> (2013)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/23054007">https://www.ncbi.nlm.nih.gov/pubmed/23054007</a>
Culture supernatant	Mouse	J.Y. Seo, <i>et al.</i> (2015)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26548345">https://www.ncbi.nlm.nih.gov/pubmed/26548345</a>
Culture supernatant	Mouse	Y.A. Bae, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27382358">https://www.ncbi.nlm.nih.gov/pubmed/27382358</a>
Culture supernatant	Mouse	C.G. Lee, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26791584">https://www.ncbi.nlm.nih.gov/pubmed/26791584</a>
Culture supernatant	Mouse	J. Lee, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26709074">https://www.ncbi.nlm.nih.gov/pubmed/26709074</a>
Culture supernatant	Mouse	C.Y. Wang, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27520377">https://www.ncbi.nlm.nih.gov/pubmed/27520377</a>
Liver tissue	Mouse	A. Moslehi, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/25183508">https://www.ncbi.nlm.nih.gov/pubmed/25183508</a>
Lung tissue	Mouse	A. Hsu, <i>et al.</i> (2007)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/17822444">https://www.ncbi.nlm.nih.gov/pubmed/17822444</a>
Ovaries	Mouse	L.E. Henkes, <i>et al.</i> (2008)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/18505843">https://www.ncbi.nlm.nih.gov/pubmed/18505843</a>
Plasma	Mouse	A. Nemeč, <i>et al.</i> (2012)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/21489584">https://www.ncbi.nlm.nih.gov/pubmed/21489584</a>
Plasma	Mouse	G.A. El Shoubaky, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26917974">https://www.ncbi.nlm.nih.gov/pubmed/26917974</a>
Serum	Mouse	O. Desy, <i>et al.</i> (2008)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/18684924">https://www.ncbi.nlm.nih.gov/pubmed/18684924</a>
Serum	Mouse	S. Imaizumi, <i>et al.</i> (2010)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/20619710">https://www.ncbi.nlm.nih.gov/pubmed/20619710</a>
Serum	Mouse	S.Y. Park, <i>et al.</i> (2010)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/20131233">https://www.ncbi.nlm.nih.gov/pubmed/20131233</a>
Serum	Mouse	Y. Zhang, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/25150882">https://www.ncbi.nlm.nih.gov/pubmed/25150882</a>
Serum	Mouse	M. Abdel-Daim, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/26796269">https://www.ncbi.nlm.nih.gov/pubmed/26796269</a>
Serum	Mouse	C.Y. Wang, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27520377">https://www.ncbi.nlm.nih.gov/pubmed/27520377</a>
Serum	Mouse	B. Shan, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28346409">https://www.ncbi.nlm.nih.gov/pubmed/28346409</a>
Serum	Mouse	G.U. Hong, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30195017">https://www.ncbi.nlm.nih.gov/pubmed/30195017</a>
Serum	Mouse	Y.H. Li, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29512688">https://www.ncbi.nlm.nih.gov/pubmed/29512688</a>
Serum	Mouse	Y. Lu, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29091299">https://www.ncbi.nlm.nih.gov/pubmed/29091299</a>
Serum	Mouse	H.A. Seong, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29700281">https://www.ncbi.nlm.nih.gov/pubmed/29700281</a>
Serum	Mouse	J. Huang, <i>et al.</i> (2019)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/31301301">https://www.ncbi.nlm.nih.gov/pubmed/31301301</a>

## TNF- $\alpha$ (rat) ELISA kit (ADI-900-086A)

For the quantitative determination of TNF- $\alpha$  in rat culture supernatants.

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Liver tissue	Rat	J. Bae, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30334699">https://www.ncbi.nlm.nih.gov/pubmed/30334699</a>
Plasma	Rat	H.C. Liu, <i>et al.</i> (2017)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/28911640">https://www.ncbi.nlm.nih.gov/pubmed/28911640</a>
Serum	Rat	S. Li, <i>et al.</i> (2014)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/24489777">https://www.ncbi.nlm.nih.gov/pubmed/24489777</a>

## VEGF (human) ELISA kit (ENZ-KIT156)

For the quantitative determination of human VEGF in serum, plasma, and culture supernatants

SAMPLE TYPE	SPECIES	REFERENCES	LINK
Culture supernatant	Human	A. Campanati, <i>et al.</i> (2016)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/27683131">https://www.ncbi.nlm.nih.gov/pubmed/27683131</a>
Culture supernatant	Human	J.S. Bok, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/30123053">https://www.ncbi.nlm.nih.gov/pubmed/30123053</a>
Culture supernatant	Human	N.G. Imir, <i>et al.</i> (2018)	<a href="https://www.ncbi.nlm.nih.gov/pubmed/29434991">https://www.ncbi.nlm.nih.gov/pubmed/29434991</a>