

AlphaLISA immunoassay kits: Limits of detection and range

AlphaLISA Limits of Detection and Range

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL201	VEGF	Vascular Endothelial Growth Factor	2.2 pg/mL	10.7 pg/mL	100,000 pg/mL	4 log
AL202	A β 1-40	Amyloid β 1-40	88 pg/mL	101 pg/mL	100,000 pg/mL	3 log
AL203	A β 1-42	Amyloid β 1-42	300 pg/mL	350 pg/mL	300,000 pg/mL	3 log
AL204	Insulin	Insulin	0.8 μ U/mL	1.3 μ U/mL	3,000 μ U/mL	3.5 log
			(mouse/rat: 239 pg/mL)			
			(bovine: 98 pg/mL)			
			(porcine: 59 pg/mL)			
AL205	IgG	Human Immunoglobulin G	0.24 ng/mL	N/A	1,000 ng/mL	3.5 log

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL206	EPO	Erythropoietin	1 mIU/mL	5.8 mIU/mL	30,000 mIU/mL	4 log
AL207	HIV p24	HIV p24	0.525 ng/mL	N/A	3,000 ng/mL	3.5 log
AL208	TNF α	Tumor Necrosis Factor alpha	2.2 pg/mL	3.6 pg/mL	30,000 pg/mL	4 log
AL209	Adiponectin	Adiponectin	45 pg/mL	32 pg/mL	1,000,000 pg/mL	4 log
AL210	CHO-P	CHO Host Cell Proteins	181 pg/mL	N/A	300,000 pg/mL	3 log
AL211	COMP	Cartilage Oligomeric Matrix Protein	1 ng/mL	1.5 ng/mL	3,000 ng/mL	3 log
AL212	EGFR	Epidermal Growth Factor receptor	127 pg/mL	76 pg/mL	1,000,000 pg/mL	3.5 log
AL213	G-CSF	Granulocyte-Colony Stimulating Factor	14 pg/mL	68 pg/mL	300,000 pg/mL	4 log
AL214	GH	Growth Hormone	5 pg/mL	13 pg/mL	100,000 pg/mL	4 log
AL215	GLP-1 (7-36 amide)	Glucagon-Like Peptide 1	27 pg/mL	65 pg/mL	30,000 pg/mL	3 log

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL216	GM-CSF	Granulocyte-Monocyte Colony Stimulating Factor	1.6 pg/mL	12 pg/mL	10,000 pg/mL	4 log
AL217	IFN- γ	Interferon gamma	11 pg/mL	8 pg/mL	100,000 pg/mL	5 log
AL218	IL10	Interleukin 10	39 pg/mL	14 pg/mL	300,000 pg/mL	4 log
AL219	IL17	Interleukin 17	14 pg/mL	20 pg/mL	300,000 pg/mL	4.5 log
AL220	IL1 β	Interleukin 1 β	0.6 pg/mL	4 pg/mL	30,000 pg/mL	4.5 log
AL221	IL2	Interleukin 2	1.8 pg/mL	22 pg/mL	100,000 pg/mL	4.5 log
AL222	IL3	Interleukin 3	11 pg/mL	13 pg/mL	300,000 pg/mL	4.5 log
AL223	IL6	Interleukin 6	1.3 pg/mL	8.6 pg/mL	30,000 pg/mL	4.5 log
AL224	IL8	Interleukin 8	1.1 pg/mL	0.9 pg/mL	30,000 pg/mL	4.5 log
AL225	Leptin	Leptin	50 pg/mL	34 pg/mL	100,000 pg/mL	3.5 log
AL226	NS0-P	NS0 Host Cell Proteins	1.6 ng/mL	N/A	1,000 ng/mL	3 log
AL227	Prolactin	Prolactin	125 pg/mL	367 pg/mL	300,000 pg/mL	3 log
AL228	PSA	Prostatic Specific Antigen	1.2 ng/mL	1.6 ng/mL	1,000 ng/mL	3 log
AL230	AFP	Alpha-fetoprotein	40.3 pg/mL	92.6 pg/mL	300,000 pg/mL	3.5 log

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL231	sAPP α	Amyloid Precursor Protein α , Secreted	61 pg/mL	N/A	1,000,000 pg/mL	4 log
AL232	sAPP β	Amyloid Precursor Protein β , Secreted	760 pg/mL	N/A	3,000,000 pg/mL	3.5 log
AL234	EPO receptor	Erythropoietin Receptor	124.8 pg/mL	135.9 pg/mL	300,000 pg/mL	3 log
AL235	ERBB2/HER2	Receptor Tyrosine-Protein Kinase ErbB-2 (HER2)	54.1 pg/mL	69.6 pg/mL	300,000 pg/mL	3.5 log
AL236	IGF1	Insulin-like Growth Factor 1	150 pg/mL	N/A	100,000 pg/mL	2.5 log
AL237	IGF2	Insulin-like Growth Factor 2	128 pg/mL (rat: 55 pg/mL)	N/A	300,000 pg/mL	3 log
AL238	IL1 α	Interleukin 1 α	1.3 pg/mL	5.9 pg/mL	30,000 pg/mL	4 log
AL239	IL12 (p70)	Interleukin 12 (p70)	4.9 pg/mL	7.4 pg/mL	300,000 pg/mL	4.5 log
AL240	IL13	Interleukin 13	3.9 pg/mL	7.8 pg/mL	100,000 pg/mL	4 log
AL241	IL18	Interleukin 18	5 pg/mL	8.9 pg/mL	300,000 pg/mL	4.5 log

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL242	MMP1	Matrix Metalloproteinase-1	82.6 pg/mL	382 pg/mL	300,000 pg/mL	3.5 log
AL243	MMP9	Matrix Metalloproteinase-9	10.9 pg/mL	49.7 pg/mL	300,000 pg/mL	4 log
AL244	CCL2/MCP1	Monocyte Chemoattractant Protein-1	3.8 pg/mL	3.5 pg/mL	100,000 pg/mL	4 log
AL245	Myeloperoxidase	Myeloperoxidase	63.8 pg/mL (mouse: 1225 pg/mL)	N/A	300,000 pg/mL	3.5 log
AL246	β -NGF	Beta-Nerve Growth Factor	32 pg/mL (mouse: 118 pg/mL)	52.3 pg/mL	300,000 pg/mL	4 log
AL249	Renin/Prorenin	Renin	7.6 pg/mL	11.2 pg/mL	300,000 pg/mL	4.5 log
AL250	tPA	Tissue Plasminogen Activator	11.1 pg/mL	21.3 pg/mL	300,000 pg/mL	4 log
AL251	VEGFB	Vascular Endothelial Growth Factor B	140 pg/mL (mouse: 233 pg/mL)	317 pg/mL	1,000,000 pg/mL	3.5 log

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL252	VEGFC	Vascular Endothelial Growth Factor C	6.2 pg/mL	18.5 pg/mL	100,000 pg/mL	4 log
AL253	VEGFD/FIGF	Vascular Endothelial Growth Factor D	29.9 pg/mL (mouse: 33 pg/mL)	78.9 pg/mL	300,000 pg/mL	4 log
AL254	sAPP α (C-term specific)	Amyloid Precursor Protein α , Soluble	76.4 pg/mL	N/A	1 000 000 pg/mL	4 log
AL255	sAPP β (high sensitivity)	Amyloid Precursor Protein β , Soluble	284 pg/mL	N/A	3 000 000 pg/mL	4 log
AL256	CA125	Mucin-16 (Ovarian carcinoma antigen CA125)	1.5 U/mL	1.3 U/mL	3 000 U/mL	3 log
AL257	CCL3 / MIP-1 α	C-C motif chemokine 3 (Macrophage inflammatory protein 1- α)	0.80 pg/mL	1.1 pg/mL	30 000 pg/mL	4.5 log

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL258	CCL4 / MIP-1 β	C-C motif chemokine 4 (Macrophage inflammatory protein 1- β)	2.8 pg/mL	2.8 pg/mL	100 000 pg/mL	4.5 log
AL259	CXCL10 / IP-10	C-X-C motif chemokine 10 (10 kDa IFN γ -induced protein)	8.6 pg/mL	13.7 pg/mL	100 000 pg/mL	4 log
AL260	CXCL11 / I-TAC	C-X-C motif chemokine 11 (Interferon inducible T-cell α chemoattractant)	1.4 pg/mL	1.2 pg/mL	30 000 pg/mL	4 log
AL261	E.coli proteins	E.coli proteins (Host cell proteins)	465 pg/mL	N/A	1 000 000 pg/mL	3 log
AL262	IgA	Immunoglobulin A	449 pg/mL	N/A	3 000 000 pg/mL	4 log
AL263	IgM	Immunoglobulin M	511 pg/mL	N/A	3 000 000 pg/mL	4 log

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL264	IFN- α	Interferon α	20.8 pg/mL	24.8 pg/mL	100 000 pg/mL	4 log
AL265	IFN- β	Interferon β	9.6 pg/mL	18.2 pg/mL	100 000 pg/mL	4 log
AL266	IL4	Interleukin 4	4 pg/mL	6.2 pg/mL	100 000 pg/mL	4 log
AL267	IL5	Interleukin 5	4.5 pg/mL	3.7 pg/mL	100 000 pg/mL	4 log
AL268	IL7	Interleukin 7	8.5 pg/mL	7.8 pg/mL	100 000 pg/mL	4 log
AL269	MMP2	Matrix Metalloproteinase-2	260 pg/mL	642 pg/mL	300 000 pg/mL	3 log
AL270	PCSK9	Proprotein convertase subtilisin/kexin 9	1056 pg/mL	2 140 pg/mL	300 000 pg/mL	2.5 log
AL272	TIMP1	Metalloproteinase inhibitor 1 (Tissue inhibitor of metalloproteinases)	520 pg/mL	255 pg/mL	100 000 pg/mL	2 log
AL273	TFF3	Trefoil factor 3	6.1 pg/mL	10.2 pg/mL	30 000 pg/mL	4 log
AL274	A β 1-15/16	Amyloid β 1-15 and 1-16	8.1 pg/mL		10 000 pg/mL	3 log

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL275	A β 1-40 (high specificity)	Amyloid β 1-40	16.1 pg/mL		30 000 pg/mL	3 log
AL276	A β 1-42 (high specificity)	Amyloid β 1-42	20.8 pg/mL		30 000 pg/mL	3 log
			(mouse/rat: 206 pg/mL)			
AL277	Alpha-2 macroglobulin	Alpha-2 macroglobulin	575 pg/mL		1 000 000 pg/mL	3 log
AL278	Caspase-3 active	active Caspase 3	126 pg/mL		300 000 pg/mL	3.5 log
AL279	CXCL1/GRO- α	C-X-C motif chemokine 1	2.4 pg/mL		30 000 pg/mL	4 log
AL280	CXCL9/MIG	C-X-C motif chemokine 9	1.6 pg/mL		100 000 pg/mL	4.5 log
AL281	HGFR/c-MET	Hepatocyte growth factor receptor	23.1 pg/mL		1 000 000 pg/mL	4.5 log
AL282	ICAM-1	Intercellular adhesion molecule 1	3.4 pg/mL		100 000 pg/mL	4 log
AL283	IL15	Interleukin 15	1.2 pg/mL		100 000 pg/mL	4.5 log

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL284	MMP3	Matrix Metalloproteinase-3	25.8 pg/mL		300 000 pg/mL	4 log
AL285	Myoglobin	Myoglobin	1 pg/mL	1.8 pg/mL	100 000 pg/mL	5 log
AL286	PAI-1	Plasminogen Activator Inhibitor-1	39 pg/mL	53 pg/mL	300 000 pg/mL	4 log
AL287	Residual Protein A	Protein A	10.8 pg/mL		30 000 pg/mL	3 log
AL288	Amyloid β 1-x	Amyloid β 1-x	4.5 pg/mL		30,000 pg/mL	4 log
AL289	CCL5/RANTES	C-C motif chemokine 5 (T-cell specific protein RANTES)	0.8 pg/mL		100 000 pg/mL	5 log
AL290	D-dimer	Fibrin D-dimer	41.3 pg/mL		1 000 000 pg/mL	4.5 log
AL291	HIV p24 (high sensitivity)	p24 (high sensitivity)	1.8 pg/mL	58.8 pg/mL	30 000 pg/mL	4.5 log
AL292	IgE	Immunoglobulin E	6.4 pg/mL	13.0 pg/mL	300 000 pg/mL	5 log
AL293	IL11	Interleukin 11	8.6 pg/mL	16.2 pg/mL	300 000 pg/mL	5 log
AL294	Serum albumin (human)	Serum albumin	104.8 pg/mL		1 000 000 pg/mL	4 log

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL295	Cardiac Troponin I	Cardiac Troponin I (c-TnI)	4.7 pg/mL	2.5 pg/mL	100 000 pg/mL	4.5 log
AL296	Amyloid β oligomers	Amyloid β oligomers	1.3 pmol/L		7 552 pmol/L	3.5 log
AL501	IFN- γ (mouse)	Interferon γ	1.1 pg/mL	1.1 pg/mL	100 000 pg/mL	5 log
AL502	IL10 (mouse)	Interleukin 10	13.5 pg/mL	12.3 pg/mL	300 000 pg/mL	4 log
AL503	IL1 β (mouse)	Interleukin 1 β	11.5 pg/mL	14.1 pg/mL	100 000 pg/mL	4 log
AL504	IL6 (mouse)	Interleukin 6	1.4 pg/mL	3.6 pg/mL	100 000 pg/mL	5 log
			(rat: 14 pg/mL)			
AL505	TNF α (mouse)	Tumor Necrosis Factor α	2 pg/mL	2.9 pg/mL	30 000 pg/mL	4 log
			(rat: 1.5 pg/mL)			
AL506	A β 1-40 (mouse/rat)	Amyloid β 1-40	31.7 pg/mL		30 000 pg/mL	3 log
AL507	A β 1-42 (mouse/rat)	Amyloid β 1-42	25.2 pg/mL		100 000 pg/mL	3.5 log
AL508	Albumin (mouse)	Albumin	716 pg/mL		1 000 000 pg/mL	3 log
AL509	CCL2/MCP-1 (mouse/rat)	C-C Motif Chemokine 2	6.9 pg/mL		100 000 pg/mL	4 log
			(rat: 4.0 pg/mL)			
AL510	C-peptide (mouse/rat)	C-peptide	11.1 pg/mL		30 000 pg/mL	3 log
			(mouse: 12 pg/mL)			

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL511	IL2 (mouse)	Interleukin 2	1.8 pg/mL		30 000 pg/mL	4 log
AL512	IL7 (mouse)	Interleukin 7	24.8 pg/mL		300 000 pg/mL	4 log
			(rat: 14 pg/mL)			
AL513	IL15 (mouse)	Interleukin 15	20.7 pg/mL		300 000 pg/mL	4 log
			(rat: 46 pg/mL)			
AL514	IL17A (mouse/rat)	Interleukin 17A	4.1 pg/mL		100 000 pg/mL	4 log
			(rat: 4.3 pg/mL)			
AL515	Adiponectin (mouse)	Adiponectin	3.1 pg/mL		100 000 pg/mL	4.5 log
AL516	CCL5/RANTES (mouse)	C-C motif chemokine 5 (T-cell specific protein RANTES)	2.6 pg/mL	3.5 pg/mL	30 000 pg/mL	4 log
AL517	GM-CSF (mouse)	Granulocyte-Macrophage Colony-Stimulating Factor	2.0 pg/mL	2.0 pg/mL	10 000 pg/mL	3.5 log
AL518	IL1 β (rat)	Interleukin 1 β	6.2 pg/mL		100 000 pg/mL	4.5 log

Catalog number	Analyte	Analyte (full name)	LDL (in buffer)	LDL (in serum)	HDL (in buffer and serum)	Dynamic Range (in buffer)
AL519	MMP9 (mouse)	Matrix Metalloproteinase-9	19.9 pg/mL	41.9 pg/mL	300 000 pg/mL	4 log
AL520	VEGFA (mouse)	Vascular Endothelial Growth Factor A	1.5 pg/mL		100 000 pg/mL	5 log

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

ASK wiki