

TECHNICAL BRIEF

AlphaLISA immunoassay kits: Species selectivity table

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL201	VEGF	Angiogenesis	human	y	x	x	c	c						
AL202	A β 1-40	Neurodegeneration	human	y	y	y	ss	ss	ss	ss	ss		ss	
AL203	A β 1-42	Neurodegeneration	human	y	y	y	ss	ss	ss	ss	ss		ss	
AL204	Insulin	Metabolism	human	y	y	y	y	y	ss	ss	ss	c	ss	
AL205	IgG	Biological	human	y	x	x						x	x	
AL206	EPO	Cancer	human	y	x	x				c				
AL207	HIV p24	Virology	N/A											
AL208	TNF α	Angiogenesis	human	y	x	x	x			c	c			
AL209	Adiponectin	Metabolism	human	y	x	x	ss	ss	ss	ss	ss			
AL210	CHO-P	Bioprocess	N/A											

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL211	COMP	Inflammation	human	y										
AL212	EGFR	Cancer	human	y										
AL213	G-CSF	Inflammation	human	y	x									
AL214	GH	Metabolism	human	y										
AL215	GLP-1 (7-36 amide)	Metabolism	human	y	y	y	y							y
AL216	GM-CSF	Inflammation	human	y	x	x			x					
AL217	IFN- γ	Inflammation	human	y	x	x			x					
AL218	IL10	Inflammation	human	y	x	x								
AL219	IL17	Inflammation	human	y	x	x								
AL220	IL1 β	Inflammation	human	y	x	x								
AL221	IL2	Inflammation	human	y	x	x								
AL222	IL3	Inflammation	human	y	x	x								
AL223	IL6	Inflammation	human	y	x	x								

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL224	IL8	Inflammation	human	y	N/A	N/A		x	x					
AL225	Leptin	Metabolism	human	y	x	x			x					
AL226	NS0-P	Bioprocess	N/A											
AL227	Prolactin	Metabolism	human	y										
AL228	PSA	Cancer	human	y										
AL230	AFP	Cancer	human	y						C				
AL231	sAPP α	Neurodegeneration	human	y			C			SS	C			
AL232	sAPP β	Neurodegeneration	human	y	SS	SS	SS		SS	SS	SS			
AL233	CRP	Inflammation	human	y	x	x			x					
AL234	EPO receptor	Cancer	human	y	x									
AL235	ERBB2 / HER2	Cancer	human	y										

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL236	IGF1	Metabolism	human	y	x	x	SS	SS	SS	SS		C	C	
AL237	IGF2	Metabolism	human	y	x	v	C	C						
AL238	IL1 α	Inflammation	human	y	x	x		x						
AL239	IL12 (p70)	Inflammation	human	y	x	x		x	x					
AL240	IL13	Inflammation	human	y	x	x				C				
AL241	IL18	Inflammation	human	y	x	x				SS				
AL242	MMP1	Cancer	human	y	x					C				
AL243	MMP9	Cancer	human	y	x					C				
AL244	CCL2 / MCP1	Inflammation	human	y	x	x			x		C			
AL245	Myeloperoxidase	Cardiovascular	human	y	y	x				C				
AL246	β -NGF	Cancer	human	y	y	x		C	C	SS	SS			

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL247	NT-proBNP	Cardiovascular	human	y										
AL248	Plasminogen	Cardiovascular	human	y	x		x			C	C			
AL249	Renin / Prorenin	Cardiovascular	human	y						SS				
AL250	tPA	Cardiovascular	human	y	x	x				C	C		x	
AL251	VEGFB	Angiogenesis	human	y	y									
AL252	VEGFC	Angiogenesis	human	y	x	x	C		C	C				
AL253	VEGFD / FIGF	Angiogenesis	human	y	y									
AL254	sAPP α (C-term specific)	Neurodegeneration	human	y	C	C	C	C	C	C	C			
AL255	sAPP β (high sensitivity)	Neurodegeneration	human	y	C	C	C	C	C	C	C			

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL256	CA125	Cancer	human	y						C				
AL257	CCL3 / MIP-1 α	Inflammation	human	y	x	x				C				
AL258	CCL4 / MIP-1 β	Inflammation	human	y	x	x				C				
AL259	CXCL10 / IP-10	Inflammation	human	y	x	x				SS				

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL260	CXCL11 / I-TAC	Cancer	human	y	x					SS				
AL261	E.coli proteins	Bioprocess	N/A											
AL262	IgA	Biological	human	y	x									
AL263	IgM	Biological	human	y	x	x	x							
AL264	IFN- α	Inflammation	human	y	x	x				C				
AL265	IFN- β	Cancer	human	y	x	x				C				
AL266	IL4	Inflammation	human	y	x	x				SS				
AL267	IL5	Inflammation	human	y	x	x				SS				
AL268	IL7	Inflammation	human	y	x	x				SS				
AL269	MMP2	Cancer	human	y	x	C	C	C	C	C				

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL270	PCSK9	Cardiovascular	human	y	x					C				
AL271	Tau	Neurodegeneration	human	y						C	C			
AL272	TIMP1	Cancer	human	y	x	x					CC			
AL273	TFF3	Cancer	human	y						C				
AL274	A β 1-15/16	Neurodegeneration	human	y			SS	SS	SS	SS				
AL275	A β 1-40 (high specificity)	Neurodegeneration	human	y	x		SS	SS	SS	SS				
AL276	A β 1-42 (high specificity)	Neurodegeneration	human	y	y		SS	SS	SS	SS				
AL277	Alpha-2 macroglobulin	Cardiovascular	human	y	x	x				C				

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL278	Caspase-3 active	Cancer	human	y	x	x				C				
AL279	CXCL1/GRO- α	Inflammation	human	y	x	x				SS				
AL280	CXCL9/MIG	Inflammation	human	y	x					SS				
AL281	HGFR/c-MET	Cancer	human	y	x					C				
AL282	ICAM-1	Cardiovascular	human	y	x	x				C				
AL283	IL15	Inflammation	human	y	x	x				SS				

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL284	MMP3	Cancer	human	y	x					C				
AL285	Myoglobin	Cardiovascular	human	y	x	x			x	C	C			
AL286	PAI-1	Cardiovascular	human	y	x	x				C				
AL287	Residual Protein A	Bioprocess	n/a											
AL288	A β 1-x	Neurodegeneration	human	y			SS	SS	SS	SS				
AL289	CCL5/RANTES	Inflammation	human	y	x	x				SS	SS			
AL290	D-dimer	Cardiovascular	human	y										
AL291	HIV p24 (High sensitivity)	Virology	n/a											
AL292	IgE	Biological	human	y										
AL293	IL11	Inflammation	human	y	y					SS				

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL294	Serum albumin (human)	Biological	human	y	x	x	x			C				
AL295	Cardiac Troponin I	Cardiovascular	human	y	x				x	SS				
AL296	A β oligomers	Neurodegeneration	human	y			SS	SS	SS	SS				
AL299	C-peptide	Metabolism	human											
AL300	FGF21	Metabolism	human											
AL301	CHO HCP (broad reactivity)	Bioprocess												
AL302	PER.C6 [®] HCP	Bioprocess		y	x	x				SS	SS			
AL303	IgG1	Biological	human	y	y		y			C	C			
AL304	IgG4	Biological	human											
AL305	p62	Autophagy	human											
AL306	LC3B	Autophagy	human	y	x	x								
AL501	IFN- γ (mouse)	Inflammation	mouse	x	y	x	x							
AL502	IL10 (mouse)	Inflammation	mouse	x	y	x								
AL503	IL1 β (mouse)	Inflammation	mouse	x	y	x								
AL504	IL6 (mouse)	Inflammation	mouse	x	y	y								
AL505	TNF α (mouse)	Inflammation	mouse	x	y	y								

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL506	A β 1-40 (mouse/rat)	Neurodegeneration	mouse/rat	x	y	y								
AL507	A β 1-42 (mouse/rat)	Neurodegeneration	mouse/rat	x	y	y								
AL508	Albumin (mouse)	Metabolism	mouse	x	y	x	x							
AL509	CCL2/MCP-1 (mouse/rat)	Inflammation	mouse/rat	x	y	y								
AL510	C-peptide (mouse/rat)	Metabolism	mouse/rat	x	y	y								
AL511	IL2 (mouse)	Inflammation	mouse	x	y	x								
AL512	IL7 (mouse)	Inflammation	mouse	x	y	y								
AL513	IL15 (mouse)	Inflammation	mouse	x	y	y								
AL514	IL17A (mouse/rat)	Inflammation	mouse/rat	x	y	y								
AL515	Adiponectin (mouse)	Metabolism	mouse	x	y	x								

Catalog number	Analyte	Therapeutic Area	Species	Human	Mouse	Rat	Bovine	Porcine	Dog	Chimpanzee	Orangutan	Goat	Rabbit	Guinea pig
AL516	CCL5/RANTES (mouse)	Inflammation	mouse	x	y	y								
AL517	GM-CSF (mouse)	Inflammation	mouse	x	y	x			x					
AL518	IL1 β (rat)	Inflammation	rat	x	x	y			x					
AL519	MMP9 (mouse)	Cancer	mouse	x	y	y								
AL520	VEGFA (mouse)	Angiogenesis	mouse		y	y		SS	x					
AL521	Leptin (mouse)	Metabolism	Mouse		y	y	x		x					