

AlphaLISA immunoassay kits: Sample matrix table

AlphaLISA Sample Matrix Table

ND: not determined

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|------------------------------------|---------|-----------------|----------------|
| AL201 | Vascular Endothelial Growth Factor | human | 3-step | 3-step |
| AL202 | Amyloid β 1-40 | human | 2-step | 2-step |
| AL203 | Amyloid β 1-42 | human | 2-step | 2-step |
| AL204 | Insulin | human | 2-step | 2-step |
| AL205 | Human Immunoglobulin G | human | 2-step | N/A |
| AL206 | Erythropoietin | human | 2-step | 2-step |
| AL207 | HIV p24 | N/A | 2-step | N/A |
| AL208 | Tumor Necrosis Factor alpha | human | 2-step | 3-step |
| AL209 | Adiponectin | human | 2-step | 2-step |
| AL210 | CHO Host Cell Proteins | N/A | 3-step | N/A |

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|--|---------|-----------------|----------------|
| AL211 | Cartilage Oligomeric Matrix Protein | human | 2-step | 2-step |
| AL212 | Epidermal Growth Factor receptor | human | 2-step | 2-step |
| AL213 | Granulocyte-Colony Stimulating Factor | human | 2-step | 2-step |
| AL214 | Growth Hormone | human | 2-step | 2-step |
| AL215 | Glucagon-Like Peptide 1 | human | 2-step | 2-step |
| AL216 | Granulocyte-Monocyte Colony Stimulating Factor | human | 3-step | 3-step |
| AL217 | Interferon gamma | human | 2-step | 3-step |
| AL218 | Interleukin 10 | human | 2-step | 3-step |
| AL219 | Interleukin 17 | human | 2-step | 2-step |
| AL220 | Interleukin 1 β | human | 2-step | 2-step |
| AL221 | Interleukin 2 | human | 2-step | 2-step |
| AL222 | Interleukin 3 | human | 2-step | 3-step |

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|--|---------|-----------------|----------------|
| AL223 | Interleukin 6 | human | 2-step | 2-step |
| AL224 | Interleukin 8 | human | 2-step | 2-step |
| AL225 | Leptin | human | 2-step | 2-step |
| AL226 | NS0 Host Cell Proteins | N/A | 3-step | N/A |
| AL227 | Prolactin | human | 2-step | 2-step |
| AL228 | Prostatic Specific Antigen | human | 2-step | 2-step |
| AL230 | Alpha-fetoprotein | human | 2-step | 2-step |
| AL231 | Amyloid Precursor Protein α , Soluble | human | 3-step | N/A |
| AL232 | Amyloid Precursor Protein β , Soluble | human | 3-step | N/A |
| AL233 | C-Reactive Protein | human | 2-step | 2-step |
| AL234 | Erythropoietin Receptor | human | 2-step | 2-step |
| AL235 | Receptor Tyrosine-Protein Kinase ErbB-2 (HER2) | human | 2-step | 2-step |

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|------------------------------------|---------|-----------------|----------------|
| AL236 | Insulin-like Growth Factor 1 | human | 2-step | N/A |
| AL237 | Insulin-like Growth Factor 2 | human | 2-step | N/A |
| AL238 | Interleukin 1 α | human | 3-step | 3-step |
| AL239 | Interleukin 12 (p70) | human | 2-step | 2-step |
| AL240 | Interleukin 13 | human | 2-step | 2-step |
| AL241 | Interleukin 18 | human | 2-step | 2-step |
| AL242 | Matrix Metalloproteinase-1 | human | 2-step | 2-step |
| AL243 | Matrix Metalloproteinase-9 | human | 2-step | 2-step |
| AL244 | Monocyte Chemoattractant Protein-1 | human | 2-step | 2-step |
| AL245 | Myeloperoxidase | human | 2-step | N/A |
| AL246 | Beta-Nerve Growth Factor | human | 2-step | 2-step |

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|--|---------|-----------------|-----------------|
| AL247 | N-terminal pro-B-type natriuretic peptide | human | 2-step | 2-step |
| AL248 | Plasminogen | human | 2-step | 2-step |
| AL249 | Renin / Prorenin | human | 2-step | 2-step |
| AL250 | Tissue Plasminogen Activator | human | 2-step | 2-step |
| AL251 | Vascular Endothelial Growth Factor B | human | 2-step | 2-step |
| AL252 | Vascular Endothelial Growth Factor C | human | 2-step | 2-step |
| AL253 | Vascular Endothelial Growth Factor D | human | 2-step | 2-step |
| AL254 | Amyloid Precursor Protein α , Soluble | human | HiConcentration | HiConcentration |
| AL255 | Amyloid Precursor Protein β , Soluble | human | HiConcentration | HiConcentration |

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|--|---------|-----------------|----------------|
| AL256 | Mucin-16 (Ovarian carcinoma antigen CA125) | human | 2-step | 2-step |
| AL257 | C-C motif chemokine 3 (Macrophage inflammatory protein 1- α) | human | 3-step | 3-step |
| AL258 | C-C motif chemokine 4 (Macrophage inflammatory protein 1- β) | human | 2-step | 2-step |
| AL259 | C-X-C motif chemokine 10 (10 kDa IFN γ -induced protein) | human | 2-step | 2-step |

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|---|---------|-----------------|-----------------|
| AL260 | C-X-C motif chemokine 11 (Interferon inducible T-cell α chemoattractant) | human | 2-step | 2-step |
| AL261 | E.coli proteins (Host cell proteins) | N/A | 3-step | 3-step |
| AL262 | Immunoglobulin A | human | 2-step | 2-step |
| AL263 | Immunoglobulin M | human | 2-step | 2-step |
| AL264 | Interferon α | human | HiConcentration | HiConcentration |
| AL265 | Interferon β | human | 3-step | 3-step |
| AL266 | Interleukin 4 | human | 3-step | 3-step |
| AL267 | Interleukin 5 | human | 2-step | 2-step |
| AL268 | Interleukin 7 | human | 2-step | 2-step |
| AL269 | Matrix Metalloproteinase-2 | human | 2-step | 2-step |

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|--|---------|-----------------|----------------|
| AL270 | Proprotein convertase subtilisin/kexin 9 | human | 2-step | 2-step |
| AL272 | Metalloproteinase inhibitor 1 (Tissue inhibitor of metalloproteinases) | human | 3-step | 3-step |
| AL273 | Trefoil factor 3 | human | 2-step | 2-step |
| AL274 | Amyloid β 1-15 and 1-16 | human | HiConc | |
| AL275 | Amyloid β 1-40 (high specificity) | human | HiConc | |
| AL276 | Amyloid β 1-12 (high specificity) | human | HiConc | |
| AL277 | Alpha-2 macroglobulin | human | 2-step | |
| AL278 | Caspase-3 active | human | HiConc | |
| AL279 | CXCL1/GRO- α | human | HiConc | |
| AL280 | CXCL9/MIG | human | HiConc | |

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|--|---------|-----------------|----------------|
| AL281 | HGFR/c-MET | human | HiConc | |
| AL282 | ICAM-1 | human | 3-step | |
| AL283 | Interleukin 15 | human | HiConc | |
| AL284 | Matrix Metalloproteinase-3 | human | HiConc | |
| AL285 | Myoglobin | human | HiConc | |
| AL286 | Plasminogen Activator Inhibitor-1 | human | HiConc | |
| AL287 | Protein A | | 3-step | |
| AL288 | Amyloid β 1-x | human | HiConc | |
| AL289 | C-C motif chemokine 5 (T-cell-specific protein RANTES) | human | HiConc | |
| AL290 | Fibrin D-dimer | human | HiConc | |
| AL291 | p24 (high sensitivity) | human | HiConc | |
| AL292 | Immunoglobulin E | human | HiConc | |
| AL293 | Interleukin 11 | human | 3-step | |
| AL294 | Serum albumin | human | HiConc | |

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|--|---------|-----------------|-------------------------|
| AL295 | Cardiac Troponin I (c-TnI) | Human | HiConc | |
| AL296 | Amyloid β oligomers | Human | HiConc | |
| AL299 | C-peptide | Human | HiConc | HiConc |
| AL300 | Fibroblast Growth Factor 21 | Human | 3-step | 3-step |
| AL301 | Chinese Hamster Ovary Cell - Host Cell Protein | | 3-step | N/A |
| AL302 | PERC6 Cell - Host Cell Protein | | 3-step | N/A |
| AL303 | Immunoglobulin G Subclass 1 | Human | 3-step | 3-step |
| AL304 | Immunoglobulin G Subclass 4 | Human | 3-step | 3-step |
| AL305 | p62, Sequestosome-1 | Human | 3-step | HiConc modified (cells) |
| AL306 | Microtubule-Associated Protein 1 Light Chain 3 Isoform B (Mammalian Homologue to Yeast Atg8) | Human | 3-step | HiConc modified (cells) |

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|--------------------------------|-----------|-----------------|----------------|
| AL501 | Interferon γ | Mouse | 2-step | 2-step |
| AL502 | Interleukin 10 | Mouse | 2-step | 2-step |
| AL503 | Interleukin 1 β | mouse | 3-step | 3-step |
| AL504 | Interleukin 6 | mouse | 3-step | 3-step |
| AL505 | Tumor Necrosis Factor α | mouse | 3-step | 3-step |
| AL506 | Amyloid β 1-40 | mouse/rat | HiConc | |
| AL507 | Amyloid β 1-42 | mouse/rat | HiConc | |
| AL508 | Albumin | mouse | 2-step | |
| AL509 | CCL2/MCP-1 | mouse/rat | 3-step | |
| AL510 | C-peptide | mouse/rat | 2-step | |
| AL511 | Interleukin 2 | mouse | 2-step | |
| AL512 | Interleukin 7 | mouse | HiConc | |
| AL513 | Interleukin 15 | mouse | HiConc | |
| AL514 | Interleukin 17A | mouse/rat | HiConc | |
| AL515 | Adiponectin | mouse | HiConc | |

| AlphaLISA Kit | Full Name | Species | Buffer Protocol | Serum Protocol |
|---------------|--|---------|-----------------|----------------|
| AL516 | C-C motif chemokine 5 (T-cell-specific protein RANTES) | mouse | 3-step | |
| AL517 | Granulocyte-Macrophage Colony-Stimulating Factor | mouse | 3-step | |
| AL518 | Interleukin 1 β | rat | 3-step | |
| AL519 | Matrix metalloproteinase-9 | mouse | 3-step | |
| AL520 | Vascular Endothelial Growth Factor A | mouse | 3-step | |
| AL521 | Leptin | mouse | HiConc | HiConc |

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