



Adipokine Related ELISAs

Factor	Cat#	Size	Species	Sample type
Adiponectin/Acrp30	ELH-Adiponectin-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
Angiopietin-1	ELH-ANGIOPIETIN1-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
Axl	ELH-AXL-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
BDNF	ELH-BDNF-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
bFGF	ELH-bFGF-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
bFGF	ELH-bFGF-001C	96 wells (12 strips x 8 wells)	Human	Cell lysate, Tissue lysate
BMP-4	ELH-BMP4-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
BMP-6	ELH-BMP6-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
BMP-7	ELH-BMP7-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
CNTF	ELH-CNTF-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
EGF	ELH-EGF-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
EGFR	ELH-EGFR-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
ENA-78	ELH-ENA78-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
E-Selectin-001	ELH-Eselectin-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
Fas	ELH-Fas-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
Fas	ELH-Fas-001C	96 wells (12 strips x 8 wells)	Human	Cell lysate, Tissue lysate
FGF-6	ELH-FGF6-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
GITR	ELH-GITR-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
GITR	ELH-GITR-001C	96 wells (12 strips x 8 wells)	Human	Cell lysate, Tissue lysate
HCC-4	ELH-HCC4-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
HGF	ELH-HGF-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
HGF	ELH-HGF-001C	96 wells (12 strips x 8 wells)	Human	Cell lysate, Tissue lysate
IFN-gamma	ELH-IFNgamma-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IFN-gamma	ELH-IFNgamma-001C	96 wells (12 strips x 8 wells)	Human	Cell lysate, Tissue lysate
IGF-BP-1	ELH-IGFBP1-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IGF-BP-2	ELH-IGFBP2-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IGF-BP-3	ELH-IGFBP3-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IGF-I	ELH-IGFI-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IL-1 beta	ELH-IL1beta-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IL-1alpha	ELH-IL1alpha-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IL-1alpha	ELH-IL1alpha-001C	96 wells (12 strips x 8 wells)	Human	Cell lysate, Tissue lysate
IL-1ra	ELH-IL1ra-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IL-1ra	ELH-IL1ra-001C	96 wells (12 strips x 8 wells)	Human	Cell lysate, Tissue lysate
IL-6	ELH-IL6-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IL-6	ELH-IL6-001C	96 wells (12 strips x 8 wells)	Human	Cell lysate, Tissue lysate
IL-8	ELH-IL8-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IL-8	ELH-IL8-001C	96 wells (12 strips x 8 wells)	Human	Cell lysate, Tissue lysate
IL-10	ELH-IL10-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IL-11	ELH-IL11-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IL-12P70	ELH-IL12P70-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
Leptin	ELH-Leptin-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
LIF	ELH-LIF-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
MCP-1	ELH-MCP1-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine

MCSF	ELH-MCSF-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
MIF	ELH-MIF-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
MIP-1-beta	ELH-MIP1beta-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
MMP-9	ELH-MMP9-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
MSP	ELH-MSP-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
PARC	ELH-PARC-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
RANTES	ELH-RANTES-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
SDF-1alpha	ELH-SDF1alpha-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
sTNF-RII	ELH-sTNFRII-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
TGFalpha	ELH-TGFalpha-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
TGF-beta-1	ELH-TGFbeta1-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
TGF-beta-2	ELH-TGFbeta2-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
TIMP-1	ELH-TIMP1-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
TIMP-2	ELH-TIMP2-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
TNF-alpha	ELH-TNFalpha-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
VCAM-1	ELH-VCAM1-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
VEGF	ELH-VEGF-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
VEGF	ELH-VEGF-001C	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
VEGF R2	ELH-VEGFR2-001	96 wells (12 strips x 8 wells)	Human	Serum, Plasma, Conditioned medium, Urine
IL-1-beta	ELM-IL1beta-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
IL-1-beta	ELM-IL1beta-001C	96 wells (12 strips x 8 wells)	Mouse	Cell lysate, Tissue lysate
IL-6	ELM-IL6-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
IL-6	ELM-IL6-001C	96 wells (12 strips x 8 wells)	Mouse	Cell lysate, Tissue lysate
IL-12p40p70	ELM-IL12p40p70-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
Leptin	ELM-Leptin-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
IFN-gamma	ELM-IFNgamma-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
IFN-gamma	ELM-IFNgamma-001C	96 wells (12 strips x 8 wells)	Mouse	Cell lysate, Tissue lysate
MCP-1	ELM-MCP1-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
SDF-1-alpha	ELM-SDF1alpha-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
TIMP-1	ELM-TIMP1-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
TIMP-1	ELM-TIMP1-001C	96 wells (12 strips x 8 wells)	Mouse	Cell lysate, Tissue lysate
TIMP-2	ELM-TIMP2-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
TNF-alpha	ELM-TNFalpha-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
TPO	ELM-TPO-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
VCAM-1	ELM-VCAM1-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
VEGF	ELM-VEGF-001	96 wells (12 strips x 8 wells)	Mouse	Serum, Plasma, Conditioned medium
VEGF	ELM-VEGF-001C	96 wells (12 strips x 8 wells)	Mouse	Cell lysate, Tissue lysate
CNTF	ELR-CNTF-001	96 wells (12 strips x 8 wells)	Rat	Serum, Plasma, Conditioned medium
IL-10	ELR-IL10-001	96 wells (12 strips x 8 wells)	Rat	Serum, Plasma, Conditioned medium
MCP-1	ELR-MCP1-001	96 wells (12 strips x 8 wells)	Rat	Serum, Plasma, Conditioned medium
MCP-1	ELR-MCP1-001C	96 wells (12 strips x 8 wells)	Rat	Cell lysate, Tissue lysate
TIMP-1	ELR-TIMP1-001	96 wells (12 strips x 8 wells)	Rat	Serum, Plasma, Conditioned medium
TIMP-1	ELR-TIMP1-001C	96 wells (12 strips x 8 wells)	Rat	Cell lysate, Tissue lysate
TNF-alpha	ELR-TNFalpha-001	96 wells (12 strips x 8 wells)	Rat	Serum, Plasma, Conditioned medium
VEGF	ELR-VEGF-001	96 wells (12 strips x 8 wells)	Rat	Serum, Plasma, Conditioned medium
VEGF	ELR-VEGF-001C	96 wells (12 strips x 8 wells)	Rat	Cell lysate, Tissue lysate