



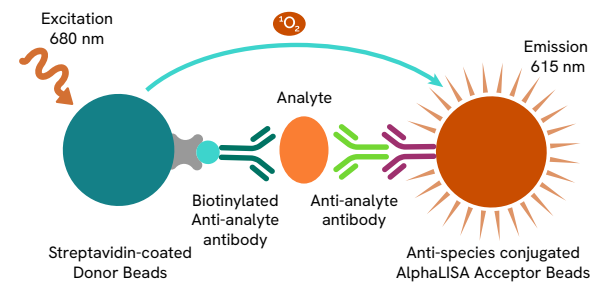
No-wash cytokine assay solutions.

Cytokines play vital roles in both innate and adaptive immunity in which they regulate immune and inflammatory responses and maintain immunological homeostasis. In addition to the immune system, they also exert diverse functions in other cells. Identification of cytokines, their receptors and signaling pathways, and application of these novel observations to the clinic has made possible the development of novel therapies for various immune diseases and related disorders.

Revvity has the tools you need to further your cytokine research with kits covering the full range of cytokines: interleukins, interferons (IFN), lymphokines, chemokines, and TNF. Our assays provide a fast alternative to ELISA or other measurement methods for researchers in the field of inflammatory disease or any other inflammation process. Benefit from our comprehensive portfolio of AlphaLISA and HTRF[®] cytokine assays that offer:

- Fast and easy homogeneous format without wash steps
- High quality results with minimal sample consumption
- Multiple cytokine measurements from a single sample
- Robust and stable assay signal for reproducible results
- Ideal for quantitative detection in cell lines like THP1 and primary cells such as PBMCs

ALPHA is a versatile, bead-based platform that enables you to assay the most complex samples in one well and with no wash steps.



HTRF is a fast, sensitive, homogeneous, and ready to use assay platform with no wash steps.

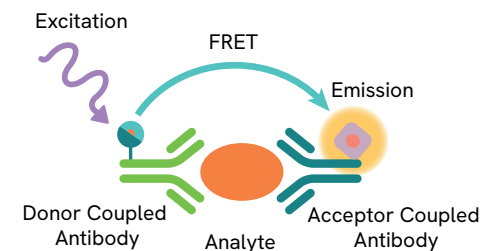


Table 1: Human cytokine assays.

	AlphaLISA	HTRF
CCL1	AL3099HV, AL3099C, AL3099F	
CCL2/MCP-1	AL244C, AL244F	62HCCL2PEG, 62HCCL2PEH
CCL3/MIP-1 α	AL257C, AL257F	62HCCL3PEG, 62HCCL3PEH
CCL4/MIP-1 β	AL258C, AL258F	62HCCL4PEG, 62HCCL4PEH
CCL5/RANTES	AL289C, AL289F	62HCCL5PEG, 62HCCL5PEH
CCL18	AL3106HV, AL3106C, AL3106F	
CCL20	62HCC20PEG, 62HCC20PEH	
CCL22	AL3188HV, AL3188C, AL3188F	62HCCL22PEG, 62HCCL22PEH
CXCL1/GRO α	AL349C, AL349F	62HCXC1PEG, 62HCXC1PEH
CXCL10	AL3163HV, AL3163C, AL3163CF	62HCX10PEG, 62HCX10PEH
CXCL11	AL260C, AL260F	
CXCL13	AL3098HV, AL3098C, AL3098F	
CXCL4/PF4	AL398HV, AL398C, AL398F	
CXCL9/MIG	AL280C, AL280F	
EPO	AL206C, AL206F	
G-CSF	AL213C, AL213F	62HGCSFPEG, 62HGCSFPEH
GM-CSF	AL216C, AL216F	62HGMCSFPEG, 62HGMCSFPEI
IFN α	AL297C, AL297F	62HIFNAPEG, 62HIFNAPEH
IFN β	AL3133C, AL3133F	62HIFNBPEG, 62HIFNBPEH
IFN γ	AL3153HV, AL3153C, AL3153F	62HIFNGPET, 62HIFNGPEG, 62HIFNGPEH
IFN λ 2	AL3031HV, AL3031C, AL3031F	
IL-1 α	AL238C, AL238F	62HIL1APEG, 62HIL1APEH
IL-1 β	AL3160HV, AL3160C, AL3160F	62HIL1BPET, 62HIL1BPEG, 62HIL1BPEH
IL-1RA	AL358HV, AL358C, AL358F	
IL-2	AL3155HV, AL3155C, AL3155F	62HIL02PET, 62HIL02PEG, 62HIL02PEH
IL-3	AL3134HV, AL3134C, AL3134F	
IL-4	AL266C, AL266F	62HIL04PEG, 62HIL04PEH
IL-5	AL267C, AL267F	62HIL05PEG, 62HIL05PEH
IL-6	AL223C, AL223F	62HIL06PET, 62HIL06PEG, 62HIL06PEH
IL-8	AL3164HV, AL3164C, AL3164F	62HIL08PET, 62HIL08PEG, 62HIL08PEH
IL-10	AL3159HV, AL3159C, AL3159F	62HIL10PEG, 62HIL10PEH
IL-11	AL293C, AL293F	62HIL11PEG, 62HIL11PEH

Table 1: Human cytokine assays continued.

	AlphaLISA	HTRF
IL-12	AL3116HV, AL3116C, AL3116F	62HIL12PEG, 62HIL12PEH
IL-12/IL-23	AL382HV, AL382C, AL382F	62HIL12P40PEG, 62HIL12P40PEH
IL-13	AL240C, AL240F	62HIL13PEG, 62HIL13PEH
IL-15	AL360HV, AL360C, AL360F	
IL-17	62HIL17PET, 62HIL17PEG, 62HIL17PEH	
IL-17A	AL3161HV, AL3161C, AL3161F	
IL-17F	AL391HV, AL391C, AL391F	
IL-18	AL3137C, AL3137F	62HIL18PEG, 62HIL18PEH
IL-21	AL3043HV, AL3043C, AL3043F	
IL-22	AL320HV, AL320C, AL320F	62HIL22PEG, 62HIL22PEH
IL-23	AL3006HV, AL3006C, AL3006F	62HIL23PEG, 62HIL23PEH
IL-25	AL3135C, AL3135F, AL3135HV	
IL-28B	AL365HV, AL365C, AL365F	
IL-31	AL3003HV, AL3003C, AL3003F	62HIL31PEG, 62HIL31PEH
IL-33	AL3042HV, AL3042C, AL3042F	
SCF	AL3019HV, AL3019C, AL3019F	
TGF β 1	AL336HV, AL336C, AL336F	
TNF α	AL3157HV, AL3157C, AL3157F	62HTNFAPET, 62HTNFAPPEG, 62HTNFAPPEH
TRAIL	AL3089HV, AL3089C, AL3089F	
TSLP	AL3040HV, AL3040C, AL3040F	
Biotin Free		
IFN γ	AL3154HV, AL3154C, AL3154F	
IL-2	AL3156HV, AL3156C, AL3156F	
IL-6	AL3025HV, AL3025C, AL3025F	
IL-8	AL3165HV, AL3165C, AL3165F	
IL-17A	AL3162HV, AL3162C, AL3162F	
TGF β 1	AL361HV, AL361C, AL361F	
TNF α	AL3158HV, AL3158C, AL3158F	

AlphaLISA available in 100 data point (HV), 500 data point (C), and 5000 data point (F) sizes.
HTRF available in 100 data point (PET), 500 data point (PEG) and 10,000 data point (PEH) sizes.

Table 2: Mouse cytokine assays.

	AlphaLISA	HTRF
CCL2/MCP1	AL509C , AL509F	62MCCL2PEG, 62MCCL2PEI
CCL4	AL3190HV, AL3190C, AL3190F	
CCL5/RANTES	AL589C , AL589F	
CXCL10/IP-10	AL556HV , AL556C , AL556F	
CXCL13	AL581HV, AL581C , AL581F	
IFN α	AL550HV , AL550C , AL550F	
IFN β	AL586HV , AL586C , AL586F	
IFN γ	AL593HV, AL593C, AL593F	62MIFNGPEG, 62MIFNGPEH
IL-1 α		62MIL1APEG, 62MIL1APEH
IL-1 β	AL592HV, AL592C, AL592F	62MIL1BPEG, 62MIL1BPEH
IL-2	AL585C , AL585F	62MIL02PEG, 62MIL02PEH
IL-3	AL587HV , AL587C , AL587F	
IL-4	AL567HV , AL567C , AL567F	
IL-5	AL569HV , AL569C , AL569F	
IL-6	AL504C , AL504F	62MIL06PEG, 62MIL06PEH
IL-10	AL502C , AL502F	
IL-13	AL565HV , AL565C , AL565F	
IL17		62MIL17PEG, 62MIL17PEH
IL-17A	AL588C , AL588F	
IL-17A homodimer	AL540HV , AL540C , AL540F	
IL-18	AL3189HV, AL3189C, AL3189F	
IL-25	AL570HV , AL570C , AL570F	
IL-33	AL566HV , AL566C , AL566F	
MCP-2	AL582HV , AL582C , AL582F	
TNF α	AL505C, AL505F	62MTNFAPEG, 62MTNFAPEH
Biotin Free		
IL-4 rhesus	AL564HV, AL564C, AL564F	
IL-6 cynomolgus, rhesus	AL555HV, AL555C, AL555F	

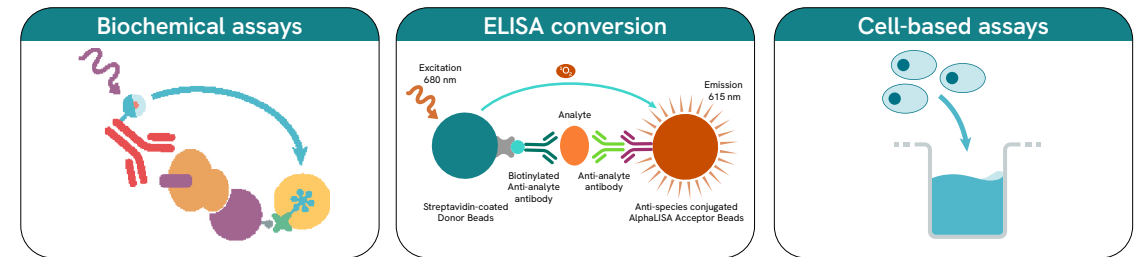
Table 3: Non-human primate cytokine assays.

	AlphaLISA	HTRF
IL-17A	AL590HV , AL590C , AL590F	
IFN γ cynomolgus	AL561HV , AL561C , AL561F	
TNF α	AL541HV , AL541C , AL541F	
TNF α cynomolgus	AL562HV , AL562C , AL562F	62CTNFAPEG, 62CTNFAPEH

AlphaLISA available in 100 data point (HV), 500 data point (C), and 5000 data point (F) sizes.
HTRF available in 100 data point (PET), 500 data point (PEG) and 10,000 data point (PEH) sizes.

Custom Labeling and Custom Assay Development Services

Don't see the assay you need listed here? Revvity partners with industry and academia professionals to develop custom solutions and shorten the drug development process using our powerful, accurate HTRF®, DELFIA® and AlphaLISA® technologies.



Contact us today: www.revvity.com

