



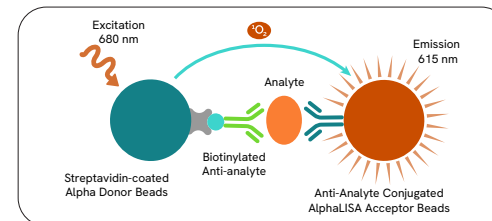
Advance your cytokine research.

The next generation of no-wash cytokine kits

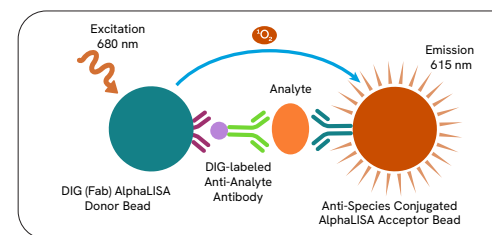
Our R&D scientists were determined to enable your research by developing new and improved versions of our AlphaLISA cytokine assays, leading to quicker, more sensitive results and faster breakthroughs. Our new AlphaLISA HIGH PERFORMANCE kits feature:

- Quantitative detection without any wash or separation steps
- Enhanced sensitivity and wide assay dynamic range
- Robust assay signal with as little as 5 μ L of sample
- Optimized performance using high quality monoclonal antibody pairs
- Results in less than 3 hours

AlphaLISA cytokine detection kits



AlphaLISA biotin free cytokine detection kits



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

New high performance kits

AlphaLISA cytokine kits	Part number
Human IFN γ	AL3153
Human IL-2	AL3155
Human TNF α	AL3157
Human IL-10	AL3159
Human IL-1 β	AL3160
Human IL-17A	AL3161
Human CXCL10	AL3163
Human IL-8	AL3164
Mouse IL-1 β	AL592
Mouse IFN γ	AL593
AlphaLISA biotin free kits	Part number
Human IFN γ	AL3154
Human IL-2	AL3156
Human TNF α	AL3158
Human IL-17A	AL3162
Human IL-8	AL3165