



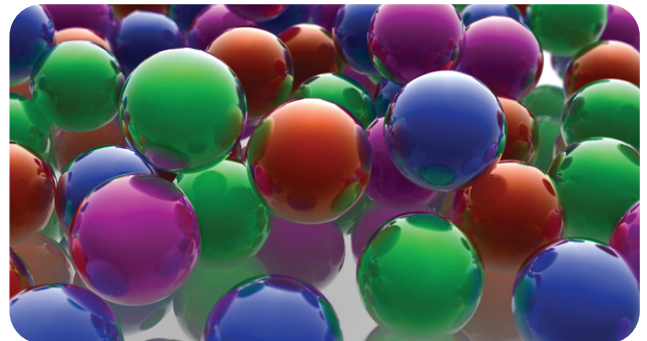
Measure more analytes in each well.

AlphaPlex technology for multiplexing assays

Now you can quantify multiple analytes in a single well for faster and more accurate results, even with low sample volumes.

AlphaPlex™ is a homogeneous, all-in-one-well multiplexing reagent technology that provides highly sensitive detection of a wide range of analytes in a simple no-wash assay format.

- High sensitivity, accuracy and well-to-well reproducibility
- Wide range of analytes: from small nucleotides and large (MDa) proteins to scarce biological samples (primary and stem cells)
- Various affinities: from pM to mM (e.g. lectin-sugar interaction)
- Cost-savings from a 5-20 times reduction in sample consumption compared to other multiplexing platforms
- Simple, no-wash assay for easier set-up, saving you time and effort
- Use with Alpha-enabled filter-based readers such as the EnVision® Multilabel Plate Reader (see document *011882_01 SHT AlphaPlex for EnVision*), and automated workstation such as JANUS®



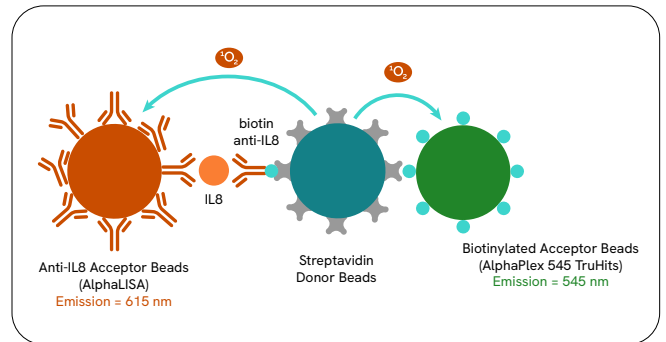
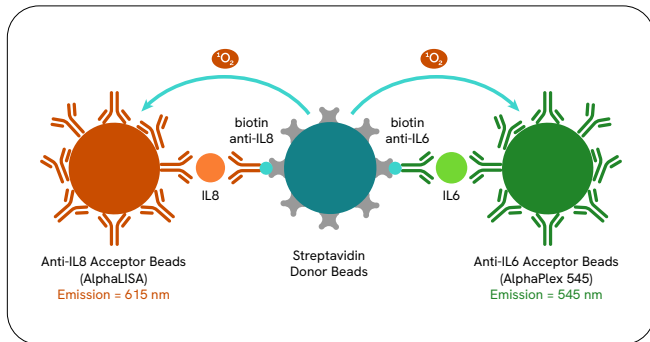
Apply the AlphaPlex reagent technology to a wide range of applications including:

- Biomarkers for PD/PK
- Biomarkers for stem cells
- Kinase: total vs phosphorylated protein
- Epigenetics markers: total histone vs specific marker
- Amyloid peptides
- IgG profiling/isotyping
- Assay normalization with housekeeping proteins

How does AlphaPlex work?

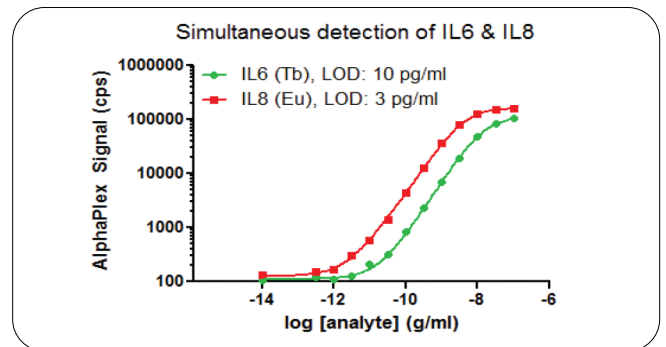
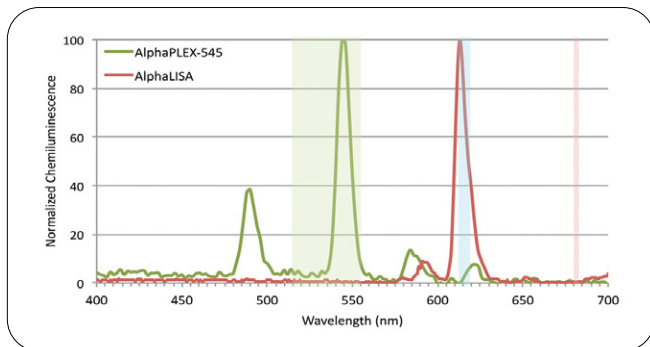
AlphaPlex is a homogeneous luminescent proximity reagent technology that allows you to quantify multiple analytes in a single well. When the Donor and Acceptor beads are brought together, a cascade of chemical reactions is set in

motion, causing a greatly amplified signal. By using multiple Acceptor beads which emit different wavelengths, multiple analytes can be detected.



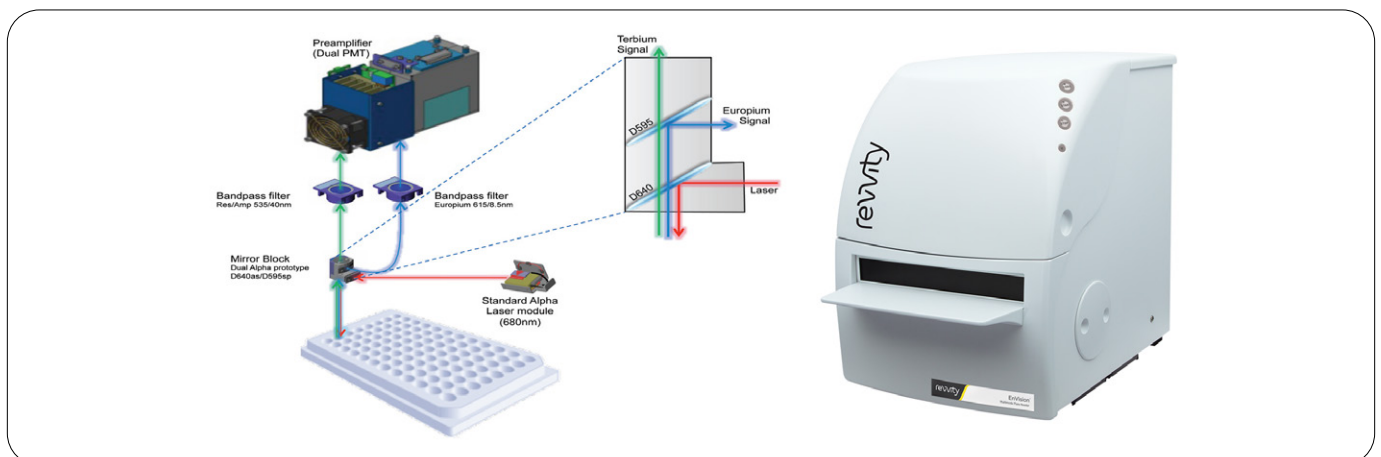
AlphaLISA IL8 and AlphaPlex 545 IL6 reagents for dual cytokine detection

AlphaLISA IL8 assay with AlphaPlex 545 TruHit control



AlphaLISA (Europium) and AlphaPlex 545 (Terbium) chemistries enable accurate quantification of analytes due to the limited spectral overlap of their respective emissions.

Biomarker pairs, such as IL6 and IL8, can be quantified with high sensitivity and selectivity using specific antibodies conjugated to AlphaPlex 545 (IL6) and AlphaLISA (IL8) Acceptor beads respectively. Sensitivity of both IL6 and IL8 assays is comparable to their best commercially available ELISA counterparts. The small error bars indicate high reproducibility/low variability.



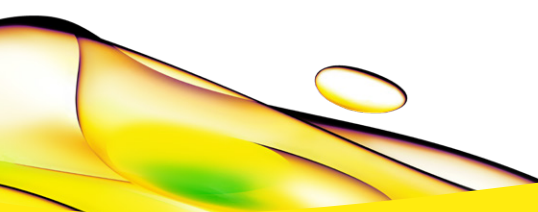
Use AlphaPlex reagent with the EnVision® Multilabel Plate Reader which is filter-based and Alpha-enabled. Please refer to document 011882_01 SHT AlphaPlex for EnVision for more details about the EnVision set-up.

Choose the product to suit your needs:

Product name	Size	Part numbers
AlphaPlex 545 Anti-Hemagglutinin Acceptor Beads	250 µg	AP170TB-C
	5 mg	AP170TB-M
	25 mg	AP170TB-R
AlphaPlex 545 Anti-Human IgG Acceptor Beads	250 µg	AP103TB-C
	5 mg	AP103TB-M
	25 mg	AP103TB-R
AlphaPlex 545 Anti-Mouse IgG Acceptor Beads	250 µg	AP105TB-C
	5 mg	AP105TB-M
	25 mg	AP105TB-R
AlphaPlex 545 Glutathione Acceptor Beads	250 µg	AP109TB-C
	5 mg	AP109TB-M
	25 mg	AP109TB-R
AlphaPlex 545 Protein A Acceptor Beads	250 µg	AP101TB-C
	5 mg	AP101TB-M
	25 mg	AP101TB-R
AlphaPlex 545 Streptavidin Acceptor Beads	250 µg	AP125TB-C
	5 mg	AP125TB-M
	25 mg	AP125TB-R
AlphaPlex 545 Unconjugated Acceptor Beads	1 mg	6782001
	5 mg	6782002
	50 mg	6782003

Product name	Size	Part numbers
AlphaPlex 545 Human IL-1β Detection Kit	100	AP220TB-HV
	500	AP220TB-C
	5,000	AP220TB-F
AlphaPlex 545 Human IL-2 Detection Kit	100	AP221TB-HV
	500	AP221TB-C
	5,000	AP221TB-F
AlphaPlex 545 Human IL-6 Detection Kit	100	AP223TB-HV
	500	AP223TB-C
	5,000	AP223TB-F
AlphaPlex 545 Human IL-17 Detection Kit	100	AP219TB-HV
	500	AP219TB-C
	5,000	AP219TB-F
AlphaPlex 545 Human and Rat IL-22 Detection Kit	100	AP320TB-HV
	500	AP320TB-C
	5,000	AP320TB-F
AlphaPlex 545 Human IFN _γ Detection Kit	100	AP217TB-HV
	500	AP217TB-C
	5,000	AP217TB-F
AlphaPlex 545 Human TNF _α Detection Kit	100	AP208TB-HV
	500	AP208TB-C
	5,000	AP208TB-F
AlphaPlex 545 TruHits Kit	500	AP900TB-D
	5,000	AP900TB-M

Contact your sales representative for more information or visit: www.revvy.com.



revvity