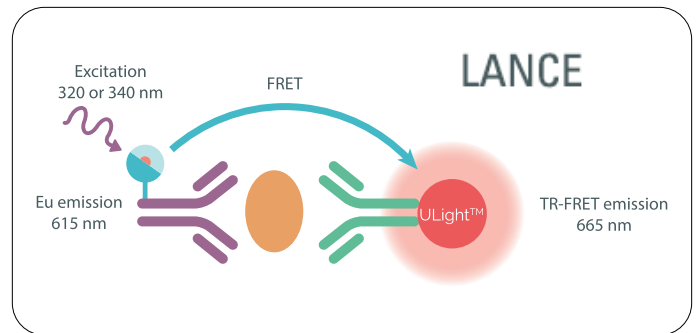
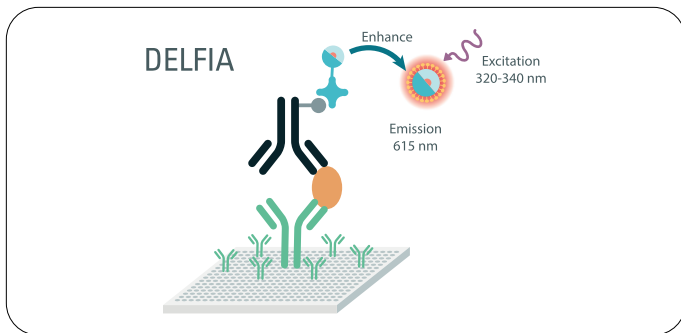
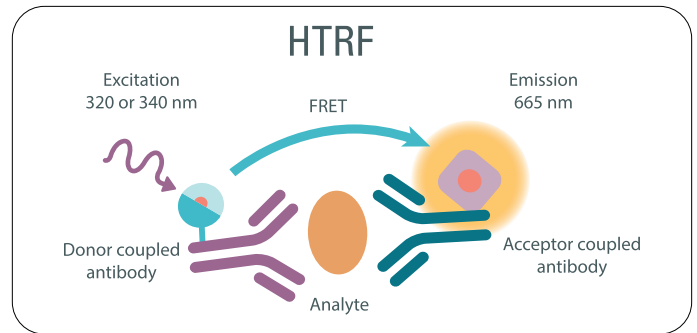
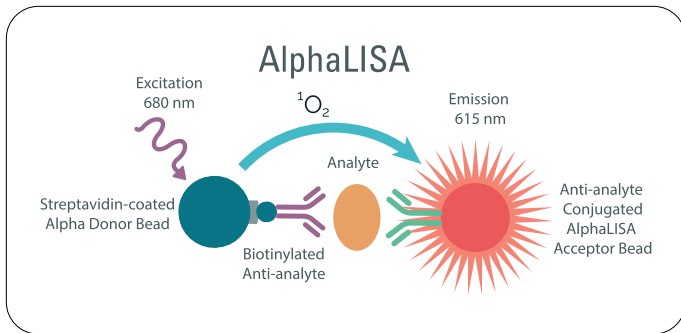
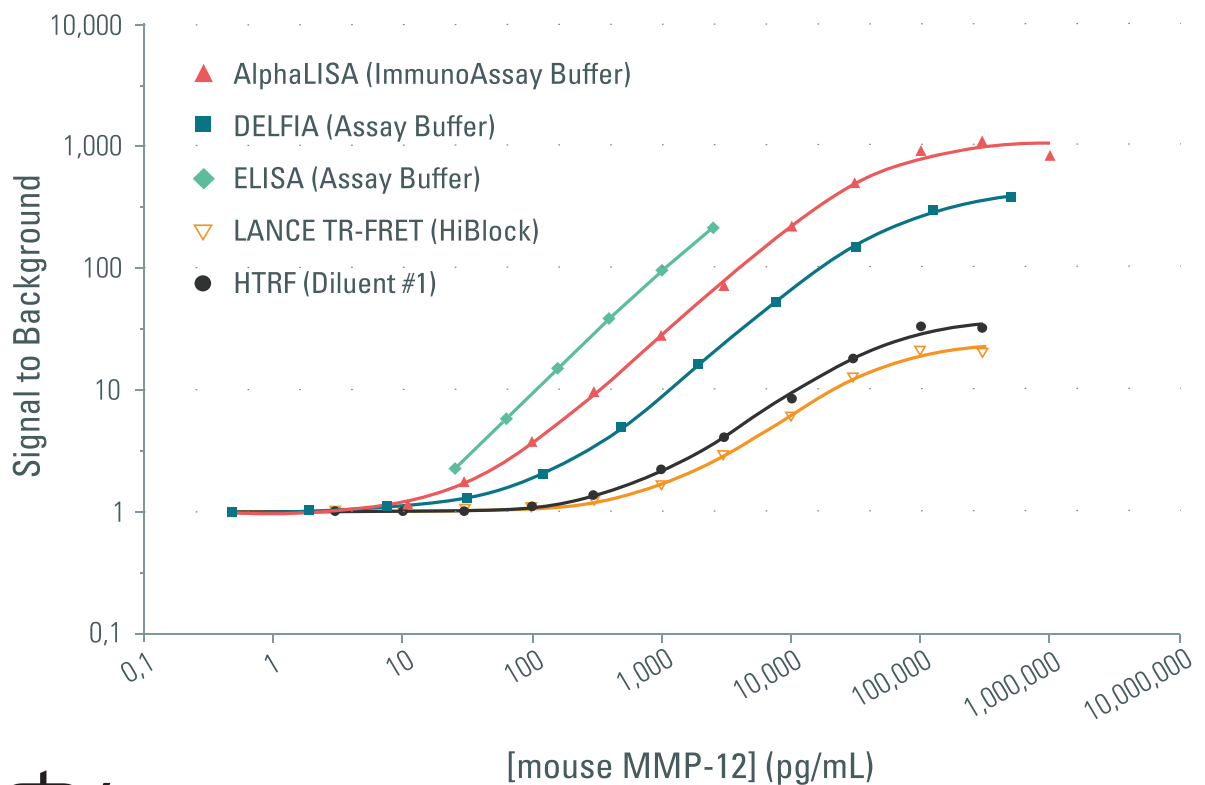


Immunoassay platforms to fit every need



Immunoassay Technologies



The right technology choice for successful assays

	ELISA	DELFI A	HTRF and LANCE Ultra	AlphaLISA
Plate Format	96-wells	96 to 384-wells	Anywhere from 96- to 1536-wells	Anywhere from 96- to 1536-wells
Miniaturization	Limited	Limited	Yes	Yes
Time To Results	4-24 hr	4-6 hr	1-2 hr	1-2 hr
			Significant time saving reallocated back to science	
Wash Steps	Yes	Yes	No	No
	While timely, in some instances wash steps can be beneficial. Wash steps may become necessary to remove nonbound analytes with higher cross-reactivity profiles and to decrease background		Absence of wash steps increases throughput, results in ease of use, reduces time to results and lowers variability	
Dynamic Range	2 Log	5 Log	4 Log	5 Log
		Large dynamic ranges are beneficial to avoid making extremely large dilutions during sample preparation. Large sample dilution are cumbersome and lead to analytical imprecision/variability		
Sensitivity	10 pg/mL	10 pg/mL	10 pg/mL	1 pg/mL
	Sufficient sensitivity to assay in serum, fluid, cell supernatant & lysates		Sufficient sensitivity to assay in cell supernatant & lysates	Sufficient sensitivity to assay in cell supernatant & lysates. May lend itself to diluting then testing serum samples, due to its superior sensitivity
Avidity	No	No	Only with streptavidin	Yes
			Avidity may be a sought after characteristic when detecting protein-protein interactions of low affinities	
Signal Stability	Up to 1 hr	Days - Years	24 hrs to days	24 hrs to days
		Allows for the plates to be left on the bench and read at a convenient time. Maximizes assay precision when incubation times vary		
Multiplexing	No	Yes (up to 3 targets)	No	Yes (up to 3 targets)
Reader Requirements	Chromogenic and Chemiluminescence Compatible	TRF Compatible	TR-FRET Compatible	Alpha Compatible
		100+ compatible reader	50+ compatible reader	10+ compatible reader
Chemistry	Enzyme-linked, HRP or AP	Fluorophore Chelate	Fluorophore Chelate (LANCE Ultra) and Fluorophore Cryptate (HTRF)	Fluorophore Embedded Beads
		Extremely robust chemistry, light & temperature insensitive	Extremely robust chemistry, light, temperature. Ion insensitive (HTRF only)	
Analyte Size	No limitation	No limitation	Subject to Förster radius (9 nm)	Subject to singlet oxygen diffusion (200 nm)
			Although LANCE, HTRF and Alpha are subject to Förster radius and singlet oxygen diffusion laws, these requirements are rarely a limitation to the development of a successful assay	

