

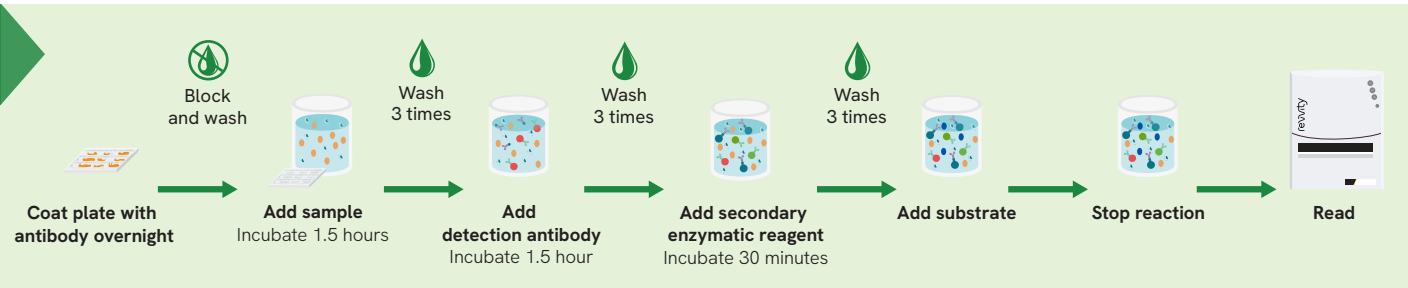
Ditch the wash steps and accelerate discovery



ELISA immunoassays are great, but what if you could improve sensitivity and increase dynamic range all while eliminating wash steps and saving time? Alpha and TR-FRET technologies do just that. These technology platforms drastically simplify the workflow of an ELISA. Just add your reagents and read - no wash needed!

ELISA

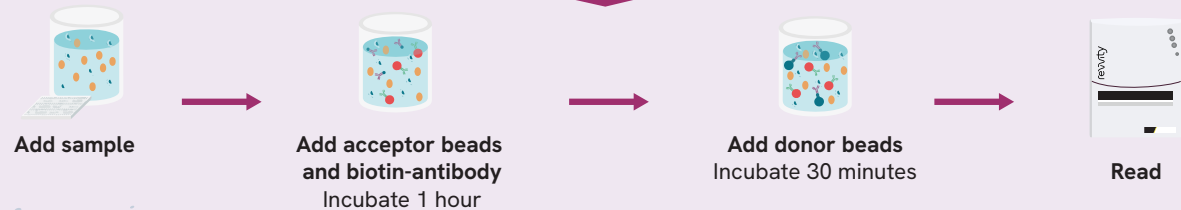
ELISA immunoassays have commonly been used for quantitation of biomarkers and other analytes in a variety of sample types.



AlphaLISA

AlphaLISA™ is a bead-based homogeneous immunoassay platform with a simple, streamlined workflow to detect and quantitate biomolecules in both simple and complex sample types.

- No washing required
- Broad sample compatibility
- Small sample size
- Wide detectable range
- Scalable



HTRF

HTRF™ is a homogeneous time-resolved fluorescence immunoassay technology that provides a simple, no wash strategy to detect and quantify proteins in as little as two hours.

- No washing required
- Saves time
- Wide detectable range
- Scalable
- Long signal stability

