

TR-FRET based Assay and Substrate.

AnaSpec, a leading provider of FRET peptides in the world, now offers custom synthesis for TR-FRET based assay development and TR-FRET substrate peptide.

- **Principle:** Use of long-lived fluorophores combined with time resolved detection (delay between excitation and emission detection) minimizes prompt fluorescence interference (this comes from proteins and compounds present in biological fluids or serum that is naturally fluorescent and is short lived).

Advantages of TR-FRET based assay.

- High sensitivity; sensitive to use at picomole level.
- Flexibility, reliability
- Use of fluorophores that has long stokes shift
- Less susceptible to interference from test compounds.
- Elimination of background fluorescence.
- Stability against photobleaching

AnaSpec: Ana Eu

Use Europium chelate, Eu (III) as donor and acceptor as QXL™ 570 or HiLiteFluor™ 647