

revvity

The easy way to approach complex biology.

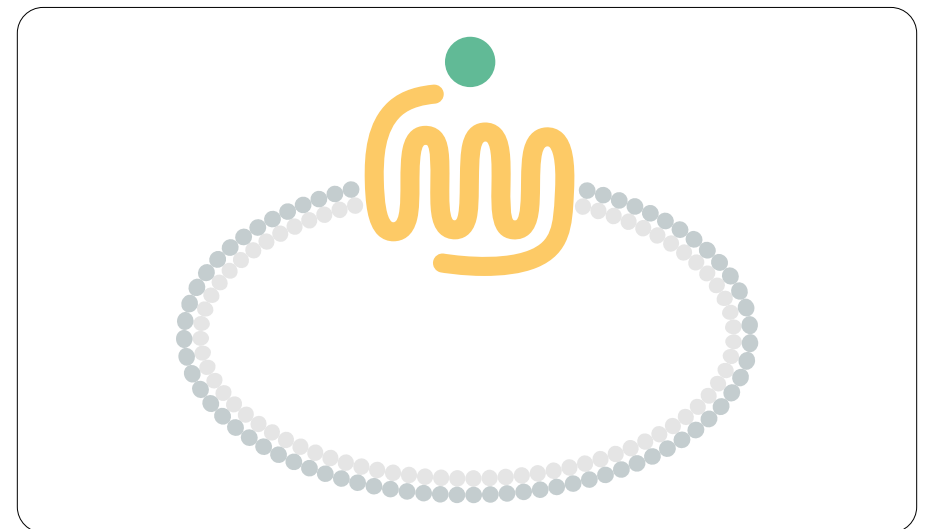
Tag-lite® binding assays

Revvity Tag-lite® binding assays offer straightforward add-and-read protocols to help you characterize the binding properties of compounds, regardless of their chemical structures (peptides, small or complex molecules), or pharmacological properties (agonist, antagonist, or inverse agonist).

Get straight to your research by running ligand binding assays directly with Revvity cells and their matching fluorescent ligands. This solution is ideal for non-radioactive, fluorescent saturation, and competitive binding assays.

Benefit from Tag-lite cellular binding assays to move away from radioactivity

- Cells and ligand pairs validated for saturation binding assays
- K_d values determination
- K_i values determination
- No effect on receptor functionality
- Frozen cells, ready to be used



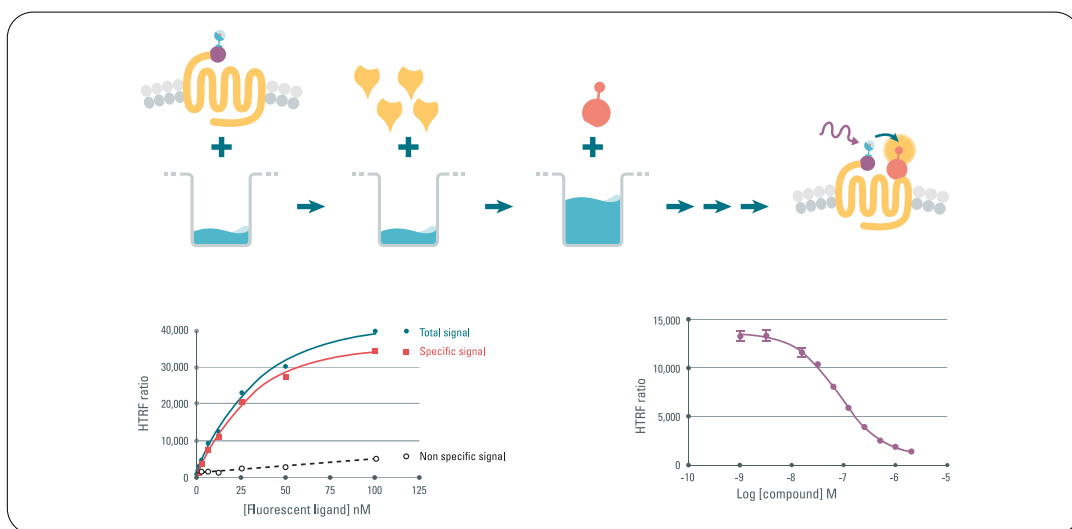
Revvity offers a wide selection of ligand binding solutions, covering over 18 GPCRs

How it works

The homogeneous alternative to radio ligand binding assays

To determine K_d , the fluorescent ligand is titrated into a solution containing a fixed amount of labeled cells, and then incubated to equilibrium. To determine K_i , the compound is titrated into a solution containing a fixed concentration of fluorescent ligand and a fixed quantity of cells. At equilibrium, the fraction of labeled ligand bound to the receptor is proportional to the recorded FRET signal. Binding affinities are calculated from this resulting signal.

Revvity provides Tag-lite transformed, and labeled cells, matching fluorescent ligands, buffer, and a complete assay procedure to guide you through assay development efficiently.



Keep your therapeutic research moving forward with faster, more robust, physiologically relevant results. Call us or visit www.revvity.com

Ordering information

| Family | Receptor | Ligand part# | Labeled cells part# | Stable cells part# |
|----------------------|-------------------|--------------|---------------------|--------------------|
| Adenosine | A1 | L0067RED | C1TT1A1 | |
| Adenosine | A2A | L0058RED | C1TT1A2A | |
| Adenosine | A2B | L0068RED | C1TT1A2B | |
| Adenosine | A3 | L0069GRE | C1TT1A3 | |
| Adrenoceptor | Beta 1 | L0023GRE | C1TT1BETA1 | |
| Adrenoceptor | Beta 2 NEW | L0011GRE | C1TT1BETA2 | C1SU1BETA2 |
| Angiotensin | AT2 | L0007RED | C1TT1AT2 | |
| Chemokine | CXCR4 NEW | L0012RED | C1TT1CXCR4 | C1SU1CXCR4 |
| Dopamine | D2 | L0002RED | C1TT1D2 | |
| Glucagon | GIPR | L0018RED | C1TT1GIPR | |
| Glucagon | GLP1 NEW | L0030RED | C1TT1GLP1 | C1SU1GLP1 |
| Orexin | OX2 | L0025RED | C1TT1OX2 | |
| Opioid | Delta | L0005RED | C1TT1DOP | |
| Opioid | Kappa | L0005RED | C1TT1KOP | |
| Opioid | Mu | L0005RED | C1TT1MOP | |
| Serotonin | 5HT1A | L0029RED | C1TT15HT1A | |
| Serotonin | 5HT4 | L0043RED | C1TT15HT4 | |
| Vasopressin/Oxytocin | V2 NEW | L0063RED | C1TT1V2 | C1PU1V2 |

Expert services to satisfy your exact needs

- In addition to providing ready-to-use solutions, Revvity's team of experienced experts also works hand-in hand with clients to create custom solutions.
- From compound characterization, to small-and large-scale reagent preparation and full assay development, Revvity is the ideal trusted partner to support all your needs.

revvity