

Tag-lite® binding assays



THE EASY WAY TO APPROACH COMPLEX BIOLOGY

Cisbio 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 Cisbio 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
- Kd values determination
- Ki values determination
- No effect on receptor functionality
- Frozen cells, ready to be used



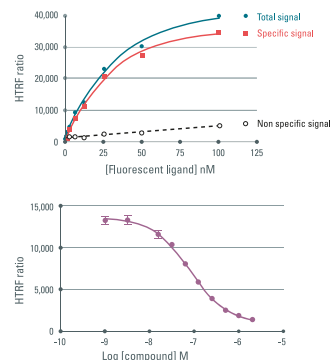
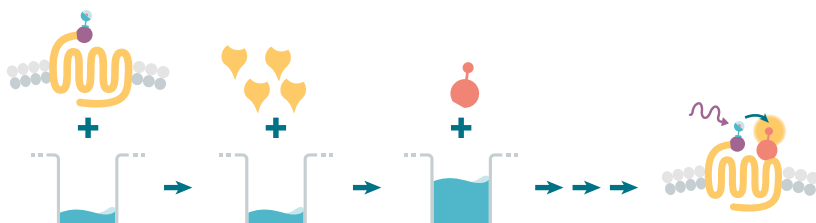
Cisbio offers a wide selection of ligand binding solutions, covering over 18 GPCRs

TAG-LITE®

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.

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



ORDERING INFORMATION

FAMILY	RECEPTOR	LIGAND PART#	LABELLED 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, Cisbio'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, Cisbio is the ideal trusted partner to support all your needs.

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

FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567

China +86 10 8080 9288 Japan +81-(0)43-306-8712

Visit www.cisbio.com to find a list of our regional distributors



cisbio