

# 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
- K<sub>d</sub> values determination
- K<sub>i</sub> values determination
- No effect on receptor functionality
- Frozen cells, ready to be used



Cisbio offers a wide selection of ligand binding solutions, covering over 42 GPCRs

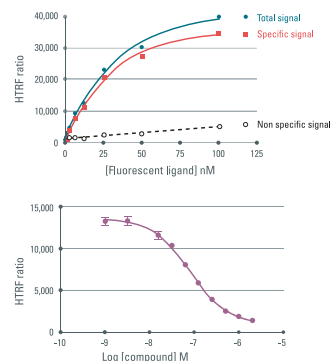
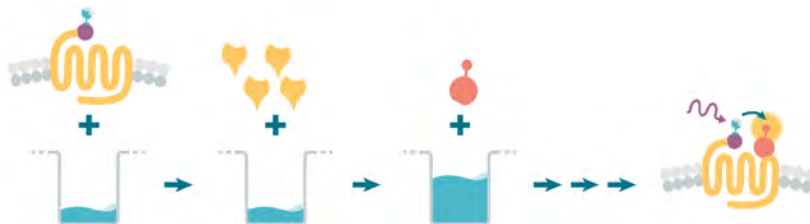
TAG-LITE®

**cisbio**  
INTERACTION IS EVERYTHING

## 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#	LABELED CELLS PART#
Adenosine	A2A	L0058RED	C1TT1A2A
Adrenoceptor	Beta 1	L0023GRE	C1TT1BETA1
Adrenoceptor	Beta 2	L0011GRE	C1TT1BETA2
Anaphylatoxin	C5L2	L0042RED	C1TT1C5L2
Angiotensin	AT2	L0007RED	C1TT1AT2
Bombesin	BB1	L0038RED	C1TT1BB1
Bombesin	BB2	L0019RED	C1TT1BB2
Bradykinin	B1	L0016RED	C1TT1B1
Bradykinin	B2	L0017RED	C1TT1B2
Chemokine	CCR5	L0024RED	C1TT1CCR5
Chemokine	CXCR4	L0012RED	C1TT1CXCR4
Cholecystokinin	CCK1	L0013RED	C1TT1CCK1
Cholecystokinin	CCK2	L0013RED	C1TT1CCK2
Class A Orphans	GPR1	L0037RED	C1TT1GPR1
Dopamine	D1	L0031GRE	C1TT1D1
Dopamine	D2	L0002RED	C1TT1D2
Dopamine	D3a	L0002RED	C1TT1D3A
Endothelin	ETA	L0022RED	C1TT1ETA
Endothelin	ETB	L0022RED	C1TT1ETB
Ghrelin	GHSR1 A	L0020RED	C1TT1GHSR1A
Glucagon	GIPR	L0018RED	C1TT1GIPR
Glucagon	GLP1	L0030RED	C1TT1GLP1

FAMILY	RECEPTOR	LIGAND PART#	LABELED CELLS PART#
Histamine	H1	L0015GRE	C1TT1H1
Histamine	H3	L0052RED	C1TT1H3
Melanocortin	MC4	L0006RED	C1TT1MC4
Melanocortin	MC5	L0006RED	C1TT1MC5
Muscarinic	M1	L0040RED	C1TT1M1
Muscarinic	M2	L0040RED	C1TT1M2
Muscarinic	M4	L0040RED	C1TT1M4
Muscarinic	M5	L0041RED	C1TT1M5
Opioid	Delta opioid	L0005RED	C1TT1DOP
Opioid	Kappa opioid	L0005RED	C1TT1KOP
Orexin	OX1	L0025RED	C1TT1OX1
Orexin	OX2	L0025RED	C1TT1OX2
Parathyroid	PTHr1	L0033RED	C1TT1PTHr1
Serotonin	5HT1A	L0029RED	C1TT15HT1A
Serotonin	5HT1B	L0045RED	C1TT15HT1B
Serotonin	5HT2C	L0066GRE	C1TT15HT2C
Serotonin	5HT4B	L0043RED	C1TT15HT4
Somatostatin	SST4	L0064RED	C1TT1SST4
Tachykinin	NK1	L0008RED	C1TT1NK1
Vasopressin/ Oxytocin	V2	L0063RED	C1TT1V2

### 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](http://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 21 5018 9880 Japan +81-(0)43-306-8712 Visit [Cisbio.com](http://Cisbio.com) to find a list of our regional distributors

