

Tag-lite Opioid receptors red antagonist - 5,000 tests

Product number: L0005RED
Batch number : 1
Storage temperature: -20°C

www.htrf.com

Product Information

For in vitro research use only

Version Mar. 11 - D

Formulation	80 µL, 10 µM solution, HEPES 50 mM pH=7.4 + 10% DMSO
Size	5,000 tests* - µ / δ / κ assays
MW	1144.4
Measure	665 nm

* 5,000 tests in a competition assay at the recommended concentration in a 384-well format (20 µL low volume)

Description

This antagonist is a naltrexone derivative labeled with a red emitting HTRF fluorescent probe.

Specificity

This ligand has been successfully used in mu-, delta- and kappa opioid Tag-lite binding assays.

Characterization

Receptor	Assay format	Kd	Working concentration*
µ	HEK293 / 384SV suspension	13 +/- 10 nM	8 nM
δ	HEK293 / 384SV suspension	5 +/- 3 nM	8 nM
κ	HEK293T / 384SV suspension	19 +/- 10 nM	8 nM

* Working concentration advised for ligand competition assays.

Storage and stability

Upon receiving the ligand, store at -20°C or colder and away from bright light.
Avoid repeated freeze-thaw cycles, as this can degrade the ligand. Individual aliquots should be prepared.

Protocole

Information available on: www.htrf.com

MSDS

Information available on: www.htrf.com