

Angiotensin receptors red ago. - 5,000 tests

Product number: L0007RED
Batch number : 1
Storage temperature: -20°C

www.htrf.com

Product Information

For in vitro research use only

Version Sept. 11 - E

Formulation	30 µL, 10 µM solution, H ₂ O + 10%MeCN
Size	5,000 tests* - AT1 / AT2 assays
MW	1670
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 agonist is an angiotensine II derivative labeled with a red emitting HTRF fluorescent probe.

Specificity

This ligand has been successfully used in AT1 and AT2 Tag-lite binding assays.

Characterization

Receptor	Assay format	Kd	Working concentration*
AT1	HEK293 / 384SV suspension	5 +/- 3 nM	3 nM
AT2	HEK293 / 384SV suspension	8 +/- 5 nM	3 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.
Centrifuge the vial before use.

Protocole

Information available on: www.htrf.com

MSDS

Information available on: www.htrf.com