

R&D, Administration and Europe Office
Cisbio Bioassays
Phone: +33 (0)4 66 79 67 05
Fax: +33 (0)4 66 79 19 20
E-mail: bioassays@cisbio.com

Japan Office
Sceti Medical Labo K.K.
Phone: +81 (0)3 5510 2932
Fax: +81 (0)3 5510 0130
E-mail: reagent@scetimedilabo.co.jp

USA Office
Cisbio US, Inc.
Phone : 888 963 4567
Fax : 781 687 1500
E-mail : htrfinfo@cisbio.us



Tag-lite Histamine H1 receptor green inv ago - 5,000 tests

Product number: L0015GRE
Batch number : 1
Storage temperature: -20°C

www.htrf.com

Product Information

For in vitro research use only
Version Oct. 10 - B

Formulation	250 µL, 20 µM solution, HEPES 50 mM pH=7.4 + 10% DMSO
Size	5,000 tests* - H1 assay
MW	874
Measure	520 nm



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

Description

This inverse agonist is a Mepyramine derivative labeled with a green emitting HTRF fluorescent probe. This product was co-developed with CellAura Technologies to be used in conjunction with Cisbio's Tag-lite® technology platform.

Specificity

This ligand has been successfully used in a Histamine H1 receptor Tag-lite binding assay.

Characterization

Receptor	Assay format	Kd	Working concentration*
H1	HEK293T / 384SV suspension	40 +/- 20nM	50 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

R&D, Administration and Europe Office

Cisbio Bioassays
Phone: +33 (0)4 66 79 67 05
Fax: +33 (0)4 66 79 19 20
E-mail: bioassays@cisbio.com

Japan Office

Sceti Medical Labo K.K.
Phone: +81 (0)3 5510 2932
Fax: +81 (0)3 5510 0130
E-mail: reagent@scetimedilabo.co.jp

USA Office

Cisbio US, Inc.
Phone : 888 963 4567
Fax : 781 687 1500
E-mail : htrfi nfo@cisbio.us



www.htrf.com

Product Information

For in vitro research use only

MSDS

Information available on: www.htrf.com