

Europe Office and Corporate Headquarters
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 Vasopressin receptors red antagonist #2 - 5,000 tests

Product number: L0001RED
Batch number : 2
Storage temperature: -20°C

www.htrf.com

Product Information

For in vitro research use only

Version Mar. 11 - D

Formulation	100 µL, 7 µM solution, H2O + 10% DMSO
Size	5,000 tests* - V2 assay
MW	1760.7
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 an Arginine-Vasopressin (AVP) derivative labeled with a red emitting HTRF fluorescent probe.

Specificity

This ligand has been successfully used both in V2R and V1AR Tag-lite binding assays.

Characterization

Receptor	Assay format	Kd	Working concentration*
V2	HEK293 / 384SV suspension	50 +/- 40 nM	5 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

Copyright © 2010 Cisbio Bioassays, France

Lumi4 is a registered trademark of Lumiphore, Inc. Lumi4-based products may be covered under one or more of U.S. patents: 6,515,113, 6,864,103, 7,442,558, 6,406,297, 7,018,850, and 7,404,912 or international counterparts.

HTRF®, Tag-lite®, and the HTRF® logo are trademarks belonging to Cisbio Bioassays.

HTRF® products are manufactured under one or more of the following patents and foreign equivalent: EP 0 180 492 / US 4,927,923 / US 5,220,012 / US 5,432,101 - EP 0 321 353 / US 5,457,185 / US 5,534,622 / US 5,346,996 / US 5,162,508 - EP 0 539 477 / US 5,512,493 - EP 0 539 435 / US 5,627,074 - EP 0 569 496 / US 5,527,684.

Cisbio Bioassays hereby grants to those buying HTRF® products from Cisbio Bioassays or its affiliates / distributors, a worldwide, non-exclusive, royalty-free, limited license to use HTRF® technology with said products for in-house life science research only. Signal amplification and correction using HTRF® technology are covered by the following U.S. patents or patent applications and foreign equivalents: US 5,512,493 - US 5,527,684 - US 6,352,672. SNAP-tag and CLIP-tag are trademarks exclusively licensed to New England Biolabs.