

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](http://www.htrf.com)

## Product Information

For in vitro research use only

Version Oct. 10 - B

## Tag-lite Beta adrenergic receptor green antagonist - 5,000 tests

**Product number:** L0011GRE  
**Batch number :** 1  
**Storage temperature:** -20°C



<b>Formulation</b>	60 µL, 10 µM solution, HEPES 50 mM pH=7.4 + 10% DMSO
<b>Size</b>	5,000 tests* - β2 assay
<b>MW</b>	760
<b>Measure</b>	520 nm

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

## Description

This beta2 adrenoceptor antagonist is a propranolol 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 Beta 2-adrenergic Tag-lite binding assay.

## Characterization

Receptor	Assay format	Kd	Working concentration*
β2	HEK293T / 384SV suspension	17 +/- 11 nM	6 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](http://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](mailto: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](mailto:reagent@scetimedilabo.co.jp)

**USA Office**

Cisbio US, Inc.  
Phone : 888 963 4567  
Fax : 781 687 1500  
E-mail : [htrfinfo@cisbio.us](mailto:htrfinfo@cisbio.us)



[www.htrf.com](http://www.htrf.com)

## Product Information

For in vitro research use only

MSDS

---

Information available on: [www.htrf.com](http://www.htrf.com)