



TAG-LITE® GLP1 RECEPTOR RED AGONIST

Product information rev 3 - March 2015

Fluorescent ligand for 5,000 tests*
Part # L0030RED**
Concentration indicated on the vial label
MW: 4999
Store at: -20°C OR BELOW

* Sufficient for 5,000 tests using a 384-well small volume plate (20 µL)

**See expiration date on product description data sheet

For research use only. Not for use in diagnostic procedures.

DESCRIPTION

This GLP1 receptor red agonist is an Exendin4 derivative labeled with a red emitting HTRF fluorescent probe.

It is suitable for both saturation binding assays (K_d) and competitive binding assays (K_i). At equilibrium, the fraction of labeled ligand bound to the receptor is proportional to the FRET signal recorded. Binding affinities can be calculated from this signal.

SPECIFICITY

This ligand has been successfully used in GLP1 Tag-lite binding assay using Glucagon GLP1 cell line Ref#C1TT1GLP1.

STORAGE AND HANDLING

Upon receiving the ligand, store at -20°C or below.

The GLP1 receptor red agonist is provided in DMSO.

NOTES:

1. We recommend centrifuging the vial of fluorescent ligand after thawing
2. Avoid repeated freeze-thaw cycles, as this can degrade the ligand. Individual aliquots should be prepared and frozen for additional use
3. We recommend the use of Tag-lite Buffer (5X concentrate), 100 mL (Cisbio Part# LABMED) for preparing working solutions
4. Specific fluorescence should be read at 665 nm
5. For more informations about Tag-lite® binding assays please refer to <http://www.cisbio.com/drug-discovery/tag-lite-binding-assays>

Copyright 2015 Cisbio Bioassays. All rights reserved. HTRF, Tag-lite, EPLigeneous and the HTRF logo are trademarks or registered trademarks of Cisbio Bioassays.

FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567

China +86-21-5018-9880 Japan +81-(0)43-306-8712

Visit www.cisbio.com to find a list of our regional distributors

