



TAG-LITE® 5HT4 RECEPTOR RED AGONIST

Product Information Revision: #03 of September 2023

Fluorescent ligand for 5,000 tests*

Part # L0043RED**

Concentration indicated on the vial label

MW: 988

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 5HT4 receptor red agonist is an ML10302 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 5HT4 Tag-lite binding assay using Serotonin 5HT4 cell line Ref#C1TT15HT4.

STORAGE AND HANDLING

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

The 5HT4 receptor red agonist is provided in 50mM Hepes buffer pH 7.4; 10% 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 (Revvity 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 www.revvity.com



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

Manufactured by Cisbio Bioassays - Parc Marcel Boiteux - 30200 Codolet - FRANCE

www.revvity.com

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.