

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT NTS1 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1276-1

**Lot Number:** C1276-1-013113

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** CHO-K1

**Transfection:** Expression vector containing full-length human NTSR1 cDNA (GenBank Accession Number NM\_002531.2) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

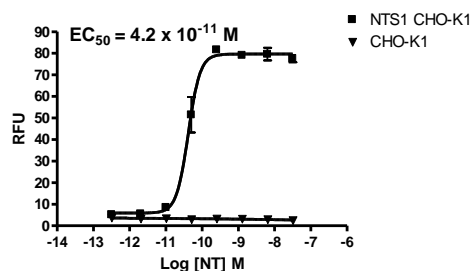
**Propagation Medium:** DMEM/F12, 10% FBS, 10  $\mu$ g/mL puromycin

**Stability:** In progress

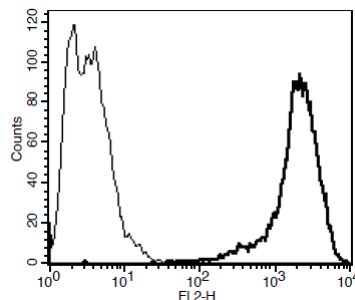
**Background:** Human NTS1 is a 418 amino acid transmembrane protein in the neurotensin receptor family. It is primarily expressed in the central nervous system and in peripheral tissues. The neurotensin family plays an important role in a wide range of biological activities, such as hypotension, hyperglycemia, and regulation of vascular permeability. Recent studies have shown that NTS1 antagonists have great potential to become novel drugs for the treatment of neurodegenerative diseases, such as Parkinson's disease.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

#### References:

- Ferraro L, et al. (2009) "Emerging evidence for neurotensin receptor 1 antagonists as novel pharmaceuticals in neurodegenerative disorders." *Mini Rev Med Chem.* 9(12):1429-38.
- Vincent JP, et al. (1999) "Neurotensin and neurotensin receptors." *Trends Pharmacol Sci.* 20(7):302-9.

Wikberg JE, et al. (2008) "Targeting melanocortin receptors: an approach to treat weight disorders and sexual disorders." *Nat Rev Drug Discov.* 7(4):307-23.

**FOR RESEARCH USE ONLY.**

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.