

PRODUCT DATASHEET

**Ready-to-Assay™ sst₃ Somatostatin
Receptor Frozen Cells****CATALOG NUMBER: HTS171RTA****CONTENTS:** Pack contains 2 vials of mycoplasma-free cells, 1 ml per vial. Fifty (50) mL of Media Component.**STORAGE:** Vials are to be stored in liquid N₂. Media Component at 4°C (-20°C for prolonged storage).**BACKGROUND**

Ready-to-Assay GPCR frozen cells are designed for simple, rapid calcium assays with no requirement for intensive cell culturing. Eurofins Discovery Services has optimized the freezing conditions to provide cells with high viability and functionality post-thaw. The user simply thaws the cells and resuspends them in media, dispenses cell suspension into assay plates and, following over night recovery, assays for calcium response.

Somatostatin is a 14 or 28 amino acid regulatory peptide that inhibits hormone secretion from the pituitary, pancreas, and other endocrine sites. A family of 6 GPCRs, sst₁, sst_{2A}, sst_{2B}, sst₃, sst₄ and sst₅, mediate the biological activity of somatostatins. The somatostatin receptors couple to G_i to inhibit cAMP production, and also increase MAP kinase signalling. Several tumors have been shown to overexpress somatostatin receptors, and binding of somatostatin to these tumor cells stimulates or inhibits proliferation, depending on the receptor subtypes expressed (Olias *et al.*, 2004). However, sst₃ appears to promote apoptosis, and expression of sst₃ was found to be lower in gastric cancer cells than in normal gastric mucosa, in proportion to susceptibility to apoptosis induced by somatostatin analogs (Sharma *et al.*, 1996; Hu *et al.*, 2004). In addition, sst₃ is expressed in smooth muscle cells and mediates gastrointestinal contractions. Cloned human sst₃-expressing cell line is made in the Chem-1 host, which supports high levels of recombinant sst₃ expression on the cell surface and contains high levels of the promiscuous G protein Gα15 to couple the receptor to the calcium signaling pathway. Thus, the cell line is an ideal tool for screening for agonists, antagonists and modulators at sst₃.

USE RESTRICTIONS

Please see User Agreement (Label License) for further details. **One such restriction is that the contents of the supplied vial(s) are limited to a single use and shall not be propagated and/or re-frozen by licensee.**

WARNINGS

For Research Use Only; Not for Use in Diagnostic Procedures
Not for Animal or Human Consumption

GMO

This product contains genetically modified organisms.
Este producto contiene organismos genéticamente modificados.
Questo prodotto contiene degli organismi geneticamente modificati.
Dieses Produkt enthält genetisch modifizierte Organismen.
Ce produit contient organismes génétiquement des modifiés.
Dit product bevat genetisch gewijzigde organismen.
Tämä tuote sisältää geneettisesti muutettuja organismeja.
Denna produkt innehåller genetiskt ändrade organismer.

APPLICATIONS

Calcium Flux Assays

APPLICATION DATA

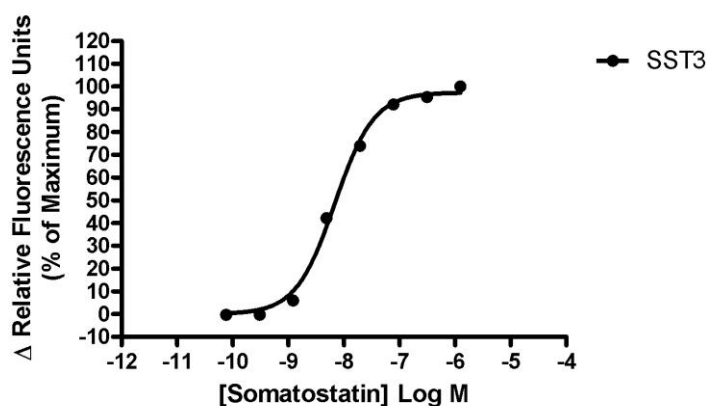


Figure 1. Representative data for activation of sst_3 receptor. Calcium flux in sst_3 -expressing Chem-1 cell line induced by Somatostatin. sst_3 -expressing Chem-1 cells were loaded with a calcium dye, and calcium flux in response to the indicated ligand(s), 4-fold serial dilution with each concentration performed in duplicate, was determined on a Molecular Devices FLIPR^{TETRA}. Maximal fluorescence signal obtained in this experiment was 4,000 RLU (Relative Light Units).

Table 1. EC₅₀ value of sst_3 -expressing Chem-1 cells.

LIGAND	ASSAY	POTENCY (nM)	REFERENCE
Somatostatin	Calcium Flux	6	Eurofins Internal Data

ASSAY SETUP

1. Immediately upon receipt, thaw cells or place cells in liquid nitrogen.
2. Thaw cells rapidly by removing from liquid nitrogen and immediately immersing in a 37°C water bath. Immediately after ice has thawed, sterilize the exterior of the vial with 70% ethanol.
3. Add 1mL of pre-warmed Media Component to each vial of cells. Place contents from two vials into a 15 mL conical tube and bring the volume to 10 mL of Media Component.
4. Centrifuge the cell suspension at 190 x g for four minutes
5. Remove supernatant and add 10.5 mL of pre-warmed Media Component to resuspend the cell pellet.
6. Seed cell suspension into appropriate assay microplate (100 μ L/well for 96-well plate, 25 μ L/well for 384-well plate).
7. When seeding is complete, place the assay plate at room temperature for 30 minutes.
8. Move assay plate to a humidified 37°C 5% CO₂ incubator for 24 hours.
9. After 24 hour incubation, remove assay plate from the incubator and wash sufficiently with Hank's Balanced Salt Solution (HBSS) supplemented with 20mM HEPES, 2.5mM Probenecid at pH 7.4 to remove

all trace of Media Component.

10. Prepare Fluo-8, AM (AAT Bioquest: 21080) Ca²⁺ dye by dissolving 1mg of Fluo-8 NW in 200 µL of DMSO. Once dissolved place 10 µL of Fluo-8 NW Ca²⁺ dye solution into 10 mL of HBSS 20mM HEPES, 2.5mM Probenecid pH 7.4 buffer and apply to assay microplate (Ca²⁺ dye at 10 µL /10 mL is sufficient for loading one (1) microplate).
11. Set-up FLIPR to dispense 3x ligand to appropriate wells in the assay plate. Set excitation wavelength at 470-495 nm (FLIPR^{TETRA}) or 485 nm (FLIPR1, FLIPR2, FLIPR3) and emission wavelength at 515-565 nm (FLIPR^{TETRA}) or emission filter for Ca²⁺ dyes (FLIPR1, FLIPR2, FLIPR3). Set pipet tip height to 5 µL below liquid level and dispense rate to 75 µL/sec (96-well format) or 50 µL/sec (384-well format). Set up plate layout and tip layout for each individual experiment. Set time course for 180 seconds, with ligand addition at 10 seconds.
12. Ligands are prepared in non-binding surface Corning plates (Corning 3605 – 96-well or Corning 3574 – 384-well).
13. After the run is complete, negative control correction is applied and data analyzed utilizing the maximum statistic.

ASSAY MATERIALS

Description	Supplier and Product Number
HBSS	Invitrogen: 14025
HEPES 1M Stock	EMD Millipore: TMS-003-C
Probenecid	Sigma: P8761
Quest Fluo-8™, AM	AAT Bioquest: 21080
Somatostatin ligand	Sigma: S9129
Non-Binding 96/384 well Plates (for ligand prep)	Corning: 3605/3574
Black (clear Bottom) cell assay plates	Corning: 3904/3712

FLIPR SETTINGS

Settings for FLIPR^{TETRA}® with ICCD camera option

Option	Setting
Read Mode	Fluorescence
Ex/Em	Ex470_495 / Em515_575
Camera Gain	2000
Gate Open	6 %
Exposure Time	0.53
Read Interval	1s
Dispense Volume	50 µl (25 µl for 384-well)
Dispense Height	25 µl (50 µl for 384-well)
Dispense Speed	75 µl L/sec (50 µl for 384-well)
Expel Volume	0 µl
Analysis	Subtract Bias Sample 1

HOST CELL

Chem-1, an adherent rat hematopoietic cell line expressing endogenous G α 15 protein

EXONGENOUS GENE EXPRESSION

SSTR3 cDNA (Accession Number: NM_001051.2; see CODING SEQUENCE below) expressed from a proprietary pHS plasmid.

CODING SEQUENCE

```

1 - ATG GAC ATG CTT CAT CCA TCA TCG GTG TCC ACG ACC TCA GAA CCT GAG AAT GCC TCC TCG GCC TGG CCC CCA - 72
1 - M D M L H P S S V S T T S E P E N A S S A W P P - 24
73 - GAT GCC ACC CTG GGC AAC GTG TCG GCA GGC CCA AGC CCG GCA GGG CTG GCC GTC AGT GGC GTT CTG ATC CCC - 144
25 - D A T L G N V S A G P S P A G L A V S G V L I P - 48
145 - CTG GTC TAC CTG GTG GTG TGC GTG GTG GGC CTG CTG GGT AAC TCG CTG GTC ATC TAT GTG GTC CTG CGG CAC - 216
49 - L V Y L V V C V V G L L G N S L V I Y V V L R H - 72
217 - ACG GCC AGC CCT TCA GTC ACC AAC GTC TAC ATC CTC AAC CTG GCG CTG GCC GAC GAG CTC TTC ATG CTG GGG - 288
73 - T A S P S V T N V Y I L N L A L A D E L F M L G - 96
289 - CTG CCC TTC CTG GCC GCC CAG AAC GCC CTG TCC TAC TGG CCC TTC GGC TCC CTC ATG TGC CGC CTG GTC ATG - 360
97 - L P F L A A Q N A L S Y W P F G S L M C R L V M - 120
361 - GCG GTG GAT GGC ATC AAC CAG TTC ACC AGC ATA TTC TGC CTG ACT GTC ATG AGC GTG GAC CGC TAC CTG GCC - 432
121 - A V D G I N Q F T S I F C L T V M S V D R Y L A - 144
433 - GTG GTA CAT CCC ACC CGC TCA GCC CGC TGG CGC ACA GCT CCG GTG GCC CGC ACG GTC AGC GCG GCT GTG TGG - 504
145 - V V H P T R S A R W R T A P V A R T V S A A V W - 168
505 - GTG GCC TCA GCC GTG GTG GTG CTG CCC GTG GTG VTC TTC TCG GGA GTG CCC CGC GGC ATG AGC ACC TGC CAC - 576
169 - L A S P A V V L P V V L P V I N L A L A D E L F M L G - 192
577 - ATG CAG TGG CCC GAG CCG GCG GCG GCC TGG CGA GCC GGC TTC ATC ATC TAC ACG GCC GCA CTG GGC TTC TTC - 648
193 - M Q W P E P A A A W R A G F I I Y T A A L G F F - 216
649 - GGG CCG CTG CTG GTC ATC TGC CTC TGC TAC CTG CTC ATC GTG GTG AAG GTG CGC TCA GCT GGG CGC CGG GTG - 720
217 - G P L C L V I C L Y L I V V K V R S A G R Y L A - 240
721 - TGG GCA CCC TCG TGC CAG CGG CGG CGG CGC TCC GAA CGC AGG GTC ACG CGC ATG GTG GTG GCG GTG GTG GCA - 792
241 - W A P S C Q R R R R S E R R V T R M V V A V V A - 264
793 - CTC TTC GTG CTC TGC TGG ATG CCC TTC TAT GTG CTC AAC ATC GTC AAC GTG GTG TGC CCA CTG CCC GAG GAG - 864
265 - L F V L C W M P F Y V L N I V N V V C P L P E E - 288
865 - CCT GCC TTC TTT GGG CTC TAC TTC CTG GTG GTG GCG CTG CCC TAT GCC AAC AGC TGT GCC AAC CCC ATC CTT - 936
289 - P A F F G L Y F L V V A L P Y A N S C A N P I L - 312
937 - TAT GGC TTC CTC TCC TAC CGC TTC AAG CAG GGC TTC CGC AGG GTC CTG CTG CGG CCC TCC CGC CGT GTG CGC -1008
313 - Y G F L S Y R F K Q G F R R V L L R P S R R V R - 336
1009 - AGC CAG GAG CCC ACT GTG GGG CCC CCG GAG AAG ACT GAG GAG GAG GAT GAG GAG GAG GAG GAT GGG GAG GAG -1080
337 - S Q E P T V G P P E K T E E E D E E E D G E E - 360
1081 - AGC AGG GAG GGG GGC AAG GGG AAG GAG ATG AAC GGC CGG GTC AGC CAG ATC ACG CAG CCT GGC ACC AGC GGG -1152
361 - S R E G G K G K E M N G R V S Q I T Q P G T S G - 384
1153 - CAG GAG CGG CCG CCC AGC AGA GTG GCC AGC AAG GAG CAG CAG CTC CTA CCC CAA GAG GCT TCC ACT GGG GAG -1224
385 - Q E R P P S R V A S K E Q Q L L P Q E A S T G E - 408
1225 - AAG TCC AGC ACG ATG CGC ATC AGC TAC CTG TGA
409 - K S S T M R I S Y L Stp

```

RELATED PRODUCTS

PRODUCT NUMBER

HTSCHEM-1RTA

HTS171M

DESCRIPTION

Ready-to-Assay™ Chem-1 host frozen cells (control cells)

ChemiScreen™ sst₃ Somatostatin receptor membrane prep

REFERENCES

1. Hu C *et al.* (2004) The effect of somatostatin and SSTR3 on proliferation and apoptosis of gastric cancer cells. *Cancer Biol Ther.* 3(8): 726-730.
2. Olias G *et al.* (2004) Regulation and function of somatostatin receptors. *J. Neurochem.* 89: 1057-1091.
3. Sharma K *et al.* (1996) Subtype-selective induction of wild-type p53 and apoptosis, but not cell cycle arrest, by human somatostatin receptor 3. *Mol. Endocrinol.* 10: 1688-1696

FOR RESEARCH USE ONLY; NOT FOR USE IN DIAGNOSTIC PROCEDURES. NOT FOR HUMAN OR ANIMAL CONSUMPTION

Unless otherwise stated in our catalog or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

No part of these works may be reproduced in any form without permission in writing.

User Agreement (Label License)

In addition to the General Terms and Conditions section, these specific terms also apply for **Ready-to-Assay™ sst₃ Somatostatin Receptor Frozen Cells, Product No. HTS171RTA**

BY USING THIS PRODUCT LICENSED TO YOU (“LICENSEE”) HEREUNDER, YOU ARE HEREBY REPRESENTING THAT YOU HAVE THE RIGHT AND AUTHORITY TO LEGALLY BIND YOURSELF OR YOUR COMPANY, AS APPLICABLE, AND ARE CONSENTING TO BE LEGALLY BOUND BY ALL OF THE TERMS OF THIS USER AGREEMENT (“AGREEMENT”). IF YOU DO NOT AGREE TO ALL THESE TERMS, DO NOT USE THE PRODUCT, AND IMMEDIATELY RETURN SUCH PRODUCTS TO THE APPLICABLE SELLER FOR A REFUND. This is a legal agreement between Licensee and Eurofins Pharma Bioanalytics Services US Inc. governing use of the Ready-to-Assay Cells products and/or any accompanying operating/use protocols (the “Product(s)”) provided to Licensee.

LICENSEE shall obtain no ownership interest in the Product or use/culture protocols accompanying the Product other than through the perpetual limited license granted herein. If the Product is licensed through an authorized Eurofins Pharma Bioanalytics Services US Inc. distributor, Licensee shall be obligated to disclose its identity to Eurofins Pharma Bioanalytics Services US Inc. to insure compliance with this User Agreement.

Limited License and Restrictions. Pursuant to the terms and conditions of this Agreement, Eurofins Pharma Bioanalytics Services US Inc. conveys to Licensee the non-exclusive and non-transferable right to use the Licensed Product only for Research Purposes conducted by Licensee (whether Licensee is an academic user or a for-profit entity). “Research Purposes” means any biological research and development application or use, including without limitation, developing, demonstrating or validating biological assays, life sciences and/or pharmaceutical research. “Research Purposes” excludes applications outside biology (including but not limited to consumer electronics or materials sciences), and specifically excludes the following applications of whatever kind or nature: Clinical Diagnostics (any use of a product or service for clinical diagnosis where data from an individual’s sample is given to such individual or used for the purpose of diagnosis or treatment of a medical condition in such individual, where that result may be used in the treatment of such individual), therapeutics, clinical imaging, environmental testing and cosmetics. Contents of the supplied vial(s) are limited to a single use and shall not be propagated and/or re-frozen by licensee. Licensee cannot sell or otherwise transfer (a) this Product or (b) materials made using this Product to a third party. Licensee may transfer information or materials made through use of this Product to a scientific collaborator, provided that such transfer is not for the commercial purposes, and that such collaborator agrees in writing: (a) not to transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for Research Purposes and not for commercial purposes. Commercial purposes means any activity by a user of the Product for consideration that may include, but is not limited to: (1) operating a service business that uses the Products to develop information or data which is resold for research and development applications; (2) use of the Product in manufacturing; (3) use of the Product for therapeutic, diagnostic or prophylactic purposes; or (4) resale of

the Product, whether or not such Product is resold for use in research. Licensee expressly represents and warrants to Eurofins Pharma Bioanalytics Services US Inc. that Licensee will properly test and use any Product purchased from Eurofins Pharma Bioanalytics Services US Inc. or its affiliated companies in accordance with the practices of a reasonable person who is an expert in the field and in strict compliance with all applicable laws and regulations, now and hereinafter enacted. Licensee agrees to comply with instructions, if any, furnished by Eurofins Pharma Bioanalytics Services US Inc. relating to the use of the Product and to not misuse the Product in any manner. Licensee shall not reverse engineer, disassemble or modify the Product or create any derivative works of the written materials accompanying the Product, including but not limited to any material data sheets or similar materials with respect to the Products' specifications. Licensee acknowledges that Eurofins Pharma Bioanalytics Services US Inc. or its affiliated companies retains ownership of all patents, copyrights, trademarks, trade secrets and other proprietary rights relating to or residing in the Product or any portion thereof.

Term and Termination. This Agreement commences upon Licensee's use of the Products, and shall remain in effect in perpetuity unless terminated sooner as set forth hereunder. Eurofins Pharma Bioanalytics Services US Inc. may terminate this Agreement immediately if Licensee breaches any provision herein. Upon any such termination, all rights granted to Licensee hereunder will immediately terminate, and Licensee shall immediately cease using the Product and, at Eurofins Pharma Bioanalytics Services US Inc.'s option, return or destroy all Products (certifying such destruction to Eurofins Pharma Bioanalytics Services US Inc. in writing).

Assignment. Licensee shall not sublicense, assign (by operation of law or otherwise) or otherwise transfer this Agreement or any of the rights or licenses granted under this Agreement without the prior written consent of Eurofins Pharma Bioanalytics Services US Inc.. Any attempted assignment, sublicense or transfer by Licensee without such consent shall be null and void.

Eurofins Pharma Bioanalytics Services US Inc. is an independent member of Eurofins Discovery Services