

## PRODUCT DATASHEET

### Ready-to-Assay™ $\alpha_{1D}$ Adrenergic Receptor Frozen Cells

#### CATALOG NUMBER: HTS216RTA

**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 N<sub>2</sub>. 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 overnight recovery, assays for calcium response.

The endogenous catecholamines epinephrine and norepinephrine have profound effects on smooth muscle activity, cardiac function, carbohydrate and fat metabolism, hormone secretion, neurotransmitter release, and central nervous system actions. These activities are mediated by GPCRs belonging to two subfamilies, the  $\alpha$ - and  $\beta$ -adrenoceptors (Bylund et al., 1994). The three members of the  $\alpha_1$  subclass of adrenoceptors,  $\alpha_{1A}$ ,  $\alpha_{1B}$  and  $\alpha_{1D}$ , couple to Gq, and promote contraction of vascular and urinary tract smooth muscle, relaxation of intestinal smooth muscle, increased contractile force in the heart, and glycogenolysis and gluconeogenesis in the liver. The different subtypes have overlapping distributions and variably contribute to these effects depending on species and tissue. The  $\alpha_{1D}$  adrenergic receptor mediates smooth muscle contraction in several tissues. In the vasculature, activation of  $\alpha_{1D}$  increases blood pressure (Tanoue et al., 2002; Hosoda et al., 2005). In the urinary tract,  $\alpha_{1D}$  promotes bladder contraction. Antagonists of  $\alpha_1$  receptors are used to treat bladder outlet obstruction, and this effect is thought to be mediated by  $\alpha_{1D}$  (Chen et al., 2005). The  $\alpha_{1D}$  adrenergic receptor has a relatively long N-terminal extracellular domain, and truncation of this domain has been shown to increase expression of the receptor at the cell surface (Pupo et al., 2003). Cloned human  $\alpha_{1D}$ -expressing cell line contains a version of  $\alpha_{1D}$  lacking residues 2-79. The cell line is made in the Chem-1 host, which supports high levels of recombinant  $\alpha_{1D}$  expression on the cell surface and contains high levels of the promiscuous G protein to couple the receptor to the calcium signaling pathway. Thus, the cell line is an ideal tool for screening for agonists and antagonists at  $\alpha_{1D}$ .

#### 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

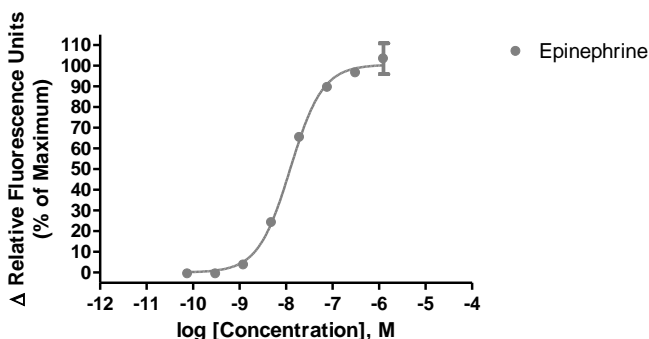


Figure 1. Representative data for activation of  $\alpha_{1D}$  receptor. Calcium flux in  $\alpha_{1D}$ -expressing Chem-1 cell line induced by Epinephrine.  $\alpha_{1D}$ -expressing Chem-1 cells were loaded with a calcium dye, and calcium flux in response to the indicated ligand(s), 4-fold serial dilution with each concentration performed in duplicate, was determined on a Molecular Devices FLIPR<sup>TETRA</sup>. Maximal fluorescence signal obtained in this experiment was 2,300 RLU (Relative Light Units).

Table 1. Comparison of EC<sub>50</sub> values of  $\alpha_{1D}$ -expressing Chem-1 cells with values described in the literature.

LIGAND	ASSAY	POTENCY (nM)	REFERENCE
Epinephrine	Calcium Flux	12	Eurofins Internal Data
Norepinephrine	Calcium Flux	3	Horie et al., 1995

### 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  $\mu$ L/well for 96-well plate, 25  $\mu$ 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% CO<sub>2</sub> 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) Ca<sup>2+</sup> dye by dissolving 1mg of Fluo-8 NW in 200  $\mu$ L of DMSO. Once dissolved place 10  $\mu$ L of Fluo-8 NW Ca<sup>2+</sup> dye solution into 10 mL of HBSS 20mM HEPES, 2.5mM Probenecid pH 7.4 buffer and apply to assay microplate (Ca<sup>2+</sup> dye at 10  $\mu$ L /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<sup>TETRA</sup>) or 485 nm (FLIPR1, FLIPR2, FLIPR3) and emission wavelength at 515-565 nm (FLIPR<sup>TETRA</sup>) or emission filter for Ca<sup>2+</sup> 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	Hyclone: SH30268.02
HEPES 1M Stock	EMD Millipore.: TMS-003-C
Probenicid	Sigma: P8761
Quest Fluo-8™, AM	AAT Bioquest: 21080
Epinephrine ligand	Sigma: E1635
Non-binding white plates (for ligand prep)	Corning: 3605(96-well)/3574(384-well)
Black (clear bottom) tissue-culture treated plates	Corning: 3904(96-well)/3712(384-well)

## FLIPR SETTINGS

Settings for FLIPR<sup>TETRA</sup>® 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 Gα15 protein.

## EXONGENOUS GENE EXPRESSION

ADRA1D cDNA (Accession Number: NM\_000678 with N-term truncation; see CODING SEQUENCE below) expressed from a proprietary PHS plasmid.

### CODING SEQUENCE

```

1 - ATG ACT TFG CGC GAT CTC CTG AGG GTC AGT TTC GAG GGA CCC CGC CGG GAC AGG AGC GCA GGG GGC TCG AGG - 72
1 - M T F R D L L S V S F E C F R P D S S A C C S S - 24
73 - CCC GCC CGC CGC CGC CGC ACC CGC CGC CGC CGC TCG GAC CGC CGC CGC CTC CGC CGC CTC CGC CGC - 144
25 - A C C C G C S A C G A A P S E C F A V G C V P G - 48
145 - GGC GCC CGC CGC CGC CGC CGC CTC CTC CGC GCA CCC AGC CGC CAC CAC AAC CGC ACC TCC CGC CGC CAG CCC - 216
49 - C A C C G C C V V G A C S C E D N R S S A C E F - 72
217 - GGG AGC GGG GGG GGG GGG GGG GAC GTG AAT GGC ACG GCG GCC GTC GGG GGA CTG GTG GTG AGC GCG CAG GGC - 288
73 - G S A G A G G D V N G T A A V G G L V V S A Q G - 96
289 - GTG GGC GTG GGC GTC TTC CTG GCA GCC TTC ATC CTT ATG GCC GTG GCA GGT AAC CTG CTT GTC ATC CTC TCA - 360
97 - V G V G V F L A A F I L M A V A G N L L V I L S - 120
361 - GTG GCC TGC AAC CGC CAC CTG CAG ACC GTC ACC AAC TAT TTC ATC GTG AAC CTG GCC GTG GCC GAC CTG CTG - 432
121 - V A C N R H L Q T V T N Y F I V N L A V A D L L - 144
433 - CTG AGC GCC ACC GTA CTG CCC TTC TCG GCC ACC ATG GAG GTT CTG GGC TTC TGG GCC TTT GGC CGC GCC TTC - 504
145 - L S A T V L P F S A T M E V L G F W A F G R A F - 168
505 - TGC GAC GTA TGG GCC GCC GTG GAC GTG CTG TGC TGC ACG GCC TCC ATC CTC AGC CTC TGC ACC ATC TCC GTG - 576
169 - C D V W A A V D V L C C T A S I L S L C T I S V - 192
577 - GAC CGG TAC GTG GGC GTG CGC CAC TCA CTC AAG TAC CCA GCC ATC ATG ACC GAG CGC AAG GGC GCC GCC ATC - 648
193 - D R Y V G V R H S L K Y P A I M T E R K A A A I - 216
649 - CTG GCC CTG CTC TGG GTC GTA GCC CTG GTG GTG TCC GTA GGG CCC CTG CTG GGC TGG AAG GAG CCC GTG CCC - 720
217 - L A L L W V V A L V V S V G P L L G W K E P V P - 240
721 - CCT GAC GAG CGC TTC TGC GGT ATC ACC GAG GAG GCG GGC TAC GCT GTC TTC TCC TCC GTG TGC TCC TTC TAC - 792
241 - P D E R F C G I T E E A G Y A V F S S V C F Y - 264
793 - CTG CCC ATG GCG GTC ATC GTG GTC ATG TAC TGC CGC GTG TAC GTG GTC GCG CGC AGC ACC ACT CGC AGC CTC - 864
265 - L P M A V I V V M Y C R V Y V V A R S T T R S L - 288
865 - GAG CCG GGC GTC AAG CGC GAG CGA GGC AAG GCC TCC GAG GTG GTG CTG CGC ATC CAC TGT CGC GGC GCG GCC - 936
289 - E A G V K R E R G K A S E V V L R I H C R G A A - 312
937 - ACG GGC GCC GAC GGG GCA CAC GGC ATG CGC AGC GCC AAG GGC CAC ACC TTC CGC AGC TCG CTC TCC GTG CGC - 1008
313 - T G G A D G A H G M R S A K G H T F R S S L S V R - 336
1009 - CTG CTC AAG TTC TCC CGT GAG AAG AAA GCG GCC AAG ACT CTG GCC ATC GTC GTG GGT GTC TTT GTG CTC TGC - 1080
337 - L L K F S R E K K A A K T L A I V V G V F V L C - 360
1081 - TGG TTC CCT TTC TTC TTT GTC CTG CCG CTC GGC TCC TTG TTC CCG CAG CTG AAG CCA TCG GAG GGC GTC TTC - 1152
361 - W F P F F F V L P L G S L F P Q L K P S E G V F - 384
1153 - AAG GTC ATC TTC TGG CTC GGC TAC TTC AAC AGC TGC GTG AAC CCG CTC ATC TAC CCC TGT TCC AGC CGC GAG - 1224
385 - K V I F W L G Y F N S C V N P L I Y P C S S R E - 408
1225 - TTC AAG CGC GCC TTC CTC CGT CTC CTG CGC TGC CAG TGC CGT CGT CGC CGG CGC CGC CCT CTC TGG CGT - 1296
409 - F K R A F L R L L R C Q C R R R R R R R P L W R - 432
1297 - GTC TAC GGC CAC CAC TGG CGG GCC TCC ACC AGC GGC CTG CGC CAG GAC TGC GCC CCG AGT TCG GGC GAC GCG - 1368
433 - V Y G H H W R A S T S G L R Q D C A P S S G D A - 456
1369 - CCC CCC GGA GCG CCG CTG GCC CTC ACC GCG CTC CCC GAC CCC GAC CCC GAA CCC CCA GGC ACG CCC GAG ATG - 1440
457 - P P G A P L A L T A L P D P D P E P P G T P E M - 480
1441 - CAG GCT CCG GTC GCC AGC CGT CGA AAG CCA CCC AGC GCC TTC CGC GAG TGG AGG CTG CTG GGG CCA TTC CGG - 1512
481 - Q A P V A S R R K P P S A F R E W R L L G P F R - 504
1513 - AGA CCC ACG ACC CAG CTG CGC GCC AAA GTC TCC AGC CTG TCG CAC AAG ATC CGC GCC GGG GGC GCG CAG CGC - 1584
505 - R P T T Q L R A K V S S L S H K I R A G G A Q R - 528
1585 - GCA GAG GCA GCG TGC GCC CAG CGC TCA GAG GTG GAG GCT GTG TCC CTA GGC GTC CCA CAC GAG GTG GCC GAG - 1656
529 - A E A A C A Q R S E V E A V S L G V P H E V A E - 552
1657 - GGC GCC ACC TGC CAG GCC TAC GAA TTG GCC GAC TAC AGC AAC CTA CGG GAG ACC GAT ATT TAA
553 - G A T C Q A Y E L A D Y S N L R E T D I Stp
    
```

## RELATED PRODUCTS

### PRODUCT NUMBER

HTSCHEM-1RTA

HTS216M

### DESCRIPTION

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

ChemiScreen™ α1D Adrenergic Family Receptor membrane prep

## REFERENCES

1. Bylund DB *et al.* (1994). IV. International Union of Pharmacology nomenclature of adrenoceptors. *Pharmacol. Rev.* 46: 121-136.
2. Chen Q *et al.* (2005) Function of the lower urinary tract in mice lacking  $\alpha_{1D}$ -adrenoceptor. *J. Urol.* 174: 370-374.
3. Horie K *et al.* (1995) Selectivity of the imidazoline  $\alpha$ -adrenoceptor agonists (oxymetazoline and cirazoline) for human cloned  $\alpha_1$ -adrenoceptor subtypes. *Br. J. Pharmacol.* 116: 1611-1618.
4. Hosoda C *et al.* (2005) Two  $\alpha_1$ -adrenergic receptor subtypes regulating the vasopressor response have differential roles in blood pressure regulation. *Mol. Pharmacol.* 67: 912-922.
5. Pupo AS *et al.* (2003) N-terminal truncation of human  $\alpha_{1D}$ -adrenoceptors increases expression of binding sites but not protein. *Eur. J. Pharmacol.* 462: 1-8.
6. Tanoue A *et al.* (2002) The  $\alpha_{1D}$ -adrenergic receptor directly regulates arterial blood pressure via vasoconstriction. *J. Clin. Invest.* 109: 765-775.

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™  $\alpha_{1D}$  Adrenergic Receptor Frozen Cells, Product No. HTS216RTA**

BY USING THE 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