

## PRODUCT DATASHEET

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

#### CATALOG NUMBER: HTS158RTA

**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. Overexpression of a constitutively active  $\alpha_{1B}$  mutant in the heart of transgenic mice resulted in cardiac hypertrophy with increased heart weight/body weight ratios. Analysis of  $\alpha_{1B}$  knock out mice has provided evidence that  $\alpha_{1B}$  is a mediator of blood pressure and aortic contractile responses induced by  $\alpha_1$  agonists (Milano et al., 1994). The locomotor and rewarding effects of psychostimulants and opiates were suppressed in mice lacking  $\alpha_{1B}$  -adrenergic receptors (Drouin et al. 2002). Cloned human  $\alpha_{1B}$ -expressing cell line is made in the Chem-1 host, which supports high levels of recombinant  $\alpha_{1B}$  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  $\alpha_{1B}$ .

#### 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

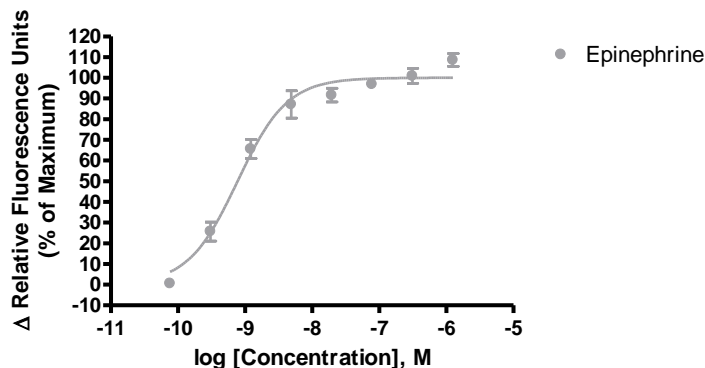


Figure 1. Representative data for activation of  $\alpha_{1B}$  receptor. Calcium flux in  $\alpha_{1B}$ -expressing Chem-1 cell line induced by epinephrine.  $\alpha_{1B}$ -expressing Chem-1 cells were loaded with a calcium dye, and calcium flux in response to the indicated ligand(s) was determined on a Molecular Devices FLIPR<sup>TETRA</sup>. Maximal fluorescence signal obtained in this experiment was 3,800 RLU (Relative Light Units).

Table 1. EC<sub>50</sub> values of  $\alpha_{1B}$ -expressing Chem-1 cells

LIGAND	ASSAY	POTENCY (nM)	REFERENCE
Epinephrine	Calcium Flux	0.8	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  $\mu$ 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 µL of DMSO. Once dissolved place 10 µL 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 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<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
Probenecid	Sigma: P8761
Quest Fluo-8™, AM	AAT Bioquest: 21080
Epinephrine ligand	Sigma: 1635
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

ADRA1B cDNA (Accession Number: NM\_000679.3; see CODING SEQUENCE below) expressed from a proprietary pHS plasmid.

## CODING SEQUENCE

```

ATG AAT CCC GAC CTG GAC ACC GGC CAC AAC ACA TCA GCA CCT GCC CAC TGG GGA GAG TTG AAA AAT GCC AAC TTC ACT
M N P D L D T G H N T S A P A H W G E L K N A N F T
GGC CCC AAC CAG ACC TCG AGC AAC TCC ACA CTG CCC CAG CTG GAC ATC ACC AGG GCC ATC TCT GTG GGC CTG GTG CTG
G P N Q T S S N S T L P Q L D I T R A I S V G L V L
GGC GCC TTC ATC CTC TTT GCC ATC GTG GGC AAC ATC CTA GTC ATC TTG TCT GTG GCC TGC AAC CGG CAC CTG CGG ACC
G A F I L F A I V G N I L V I L S V A C N R H L R T
CCC ACC AAC TAC TTC ATT GTC AAC CTG GCC ATG GCC GAC CTG CTG TTG AGC TTC ACC GTC CTG CCC TTC TCA GCG GCC
P T N Y F I V N L A M A D L L L S F T V L P F S A A
CTA GAG GTG CTC GGC TAC TGG GTG CTG GGG CGG ATC TTC TGT GAC ATC TGG GCA GCC GTG GAT GTC CTG TGC TGC ACA
L E V L G Y W V L G R I F C D I W A A V D V L C C T
GCG TCC ATT CTG AGC CTG TGC GCC ATC TCC ATC GAT CGC TAC ATC GGG GTG CGC TAC TCT CTG CAG TAT CCC ACG CTG
A S I L S L C A I S I D R Y I G V R Y S L Q Y P T L
GTC ACC CGG AGG AAG GCC ATC TTG GCG CTG CTC AGT GTC TGG GTC TTG TCC ACC GTC ATC TCC ATC GGG CCT CTC CTT
V T R I V R K A I L A L S V W V L S T V I S I G P L L
GGG TGG AAG GAG CCG GCA CCC AAC GAT GAC AAG GAG TGC GGG GTC ACC GAA GAA CCC TTC TAT GCC CTC TTC TCC TCT
G W K E P A P N D D K E C G V T E E P F Y A L F S S
CTG GGC TCC TTC TAC ATC CCT CTG GCG GTC ATT CTA GTC ATG TAC TGC CGT GTC TAT ATA GTG GCC AAG AGA ACC ACC
L G S F Y I P L A V I L V M Y C R V Y I V A K R T T
AAG AAC CTA GAG GCA GGA GTC ATG AAG GAG ATG TCC AAC TCC AAG GAG CTG ACC CTG AGG ATC CAT TCC AAG AAC TTT
K N L E A G V M K E M S N S K E L T L R I H S K N F
CAC GAG GAC ACC CTT AGC AGT ACC AAG GCC AAG GGC CAC AAC CCC AGG AGT TCC ATA GCT GTC AAA CTT TTT AAG TTC
H E D T L S S T K A K G H N P R S S I A V K L F K F
TCC AGG GAA AAG AAA GCA GCT AAG ACG TTG GGC ATT GTG GTC GGT ATG TTC ATC TTG TGC TGG CTA CCC TTC TTC ATC
S R E K K A A K T L G I V V G M F I L C W L P F F I
GCT CTA CCG CTT GGC TCC TTG TTC ACC CTG AAG CCC CCC GAC GCC GTG TTC AAG GTG GTG TTC TGG CTG GGC TAC
A L P L G S L F S T L K P P D A V F K V V F W L G Y
TTC AAC AGC TGC CTC AAC CCC ATC ATC TAC CCA TGC TCC AGC AAG GAG TTC AAG CGC GCT TTC GTG CGC ATC CTC GGG
F N S C L N P I I Y P C S S K E F K R A F V R I L G
TGC CAG TGC CGC GGC CGC GGC CGC CGA CGC CGC CGC CGT CGC CTG GGC GGC TGC GCC TAC ACC TAC CGG CCG
C Q C R G R G R R R R R R R R R R R R L G G C A Y T Y R P
TGG ACG CGC GGC GGC TCG CTG GAG CGC TCG CAG TCG CGC AAG GAC TCG CTG GAC GAC AGC GGC AGC TGC CTG AGC GGC
W T R G G S L E R S Q S R K D S L D D S G S C L S G
AGC CAG CGG ACC CTG CCC TCG GCC TCG CCG AGC CCG GGC TAC CTG GGC CGC GGC CCA CCG CCA GTC GAG CTG TGC
S Q R R T L P S A S P S P G Y L G R G A P P P V E L C
GCC TTC CCC GAG TGG AAG GCG CCC GGC GCC CTC CTG AGC CTG CCC GCG CCT GAG CCC CCC GGC CGC GGC CGC CAC
A F P E W K A P G A L L S L P A P E P P G R R G R H
GAC TCG GGC CCG CTC TTC ACC TTC AAG CTC CTG ACC GAG CCC GAG AGC CCC GGG ACC GAC GGC GGC GCC AGC AAC GGA
D S G P L F T F K L L T E P E S P G T D G G A S N G
GGC TGC GAG GCC GCG GCC GAC GTG GCC AAC GGG CAG CCG GGC TTC AAA AGC AAC ATG CCC CTG GCG CCC GGG CAG TTT
G C E A A A D V A N G Q P G F K S N M P L A P G Q F
TGA
Stop

```

## RELATED PRODUCTS

### PRODUCT NUMBER

### DESCRIPTION

**HTSCHEM-1RTA**

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

**HTS158M**

ChemiScreen™ α<sub>1B</sub> 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. Cavalli A *et al.* (1997) Decreased blood pressure response in mice deficient of the  $\alpha_{1B}$ -AR. *Proc. Natl. Acad. Sci. USA* 94: 11589–11594.
3. Milano CA *et al.* (1994) Myocardial expression of a constitutively active  $\alpha_{1B}$ -adrenergic receptor in transgenic mice induces cardiac hypertrophy. *Proc. Natl. Acad. Sci. USA* 91: 10109-10113.

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™  $\alpha_{1B}$  Adrenergic Family Receptor Frozen Cells, Product No. HTS158RTA**

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