

PRODUCT DATASHEET

Ready-to-Assay™ IP1
Prostanoid Receptor Frozen Cells

CATALOG NUMBER: HTS131RTA

Lot: 01022016

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. The freezing conditions have been optimized 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.

Prostacyclin (PGI₂) is released by vascular endothelial cells and serves as a potent vasodilator, inhibitor of platelet aggregation, and moderator of vascular smooth muscle cell proliferation–migration–differentiation (Narumiya et al. 1999). The function of prostacyclin is mediated via a seven transmembrane GPCR, IP1, which is known to couple to G_s and G_q signaling pathways. Mice lacking the IP1 receptor have shown increased susceptibility to thrombosis (Murata et al. 1997), enhanced injury-induced vascular proliferation and platelet activation (Cheng et al. 2002), as well as reperfusion injury (Xiao et al. 2001). The recent world-wide withdrawal of selective COX-2 inhibitors, rofecoxib (Vioxx™) and valdecoxib (Bextra™), is also due to their discriminating suppression of COX-2-derived prostacyclin and IP1-mediated cardioprotective effects, leading to increased risk of cardiovascular events (Fitzgerald 2004). The cloned human IP1-expressing cell line is made in the Chem-1 host, which supports high levels of recombinant IP1 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 antagonists of interactions between IP1 and its ligands.

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

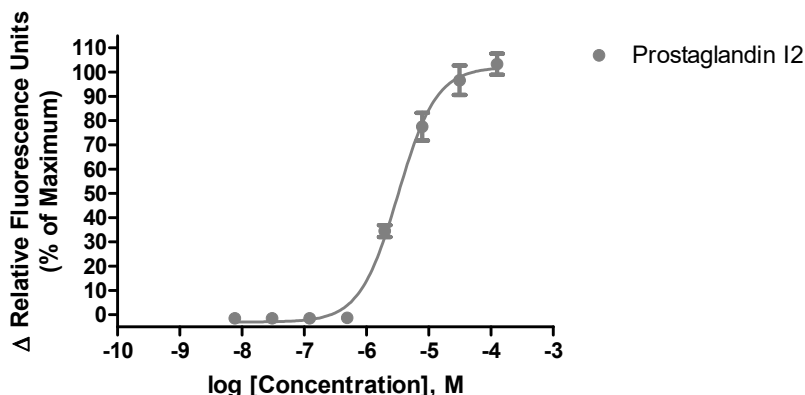


Figure 1. Representative data for activation of IP1 receptor. Calcium flux in IP1-expressing Chem-1 cell line induced by Prostaglandin I2. IP1-expressing Chem-1 cells were loaded with a calcium dye, and calcium flux in response to the indicated ligand(s), 4-fold serial dilution with each concentration performed in duplicate, was determined on a Molecular Devices FLIPR^{TETRA} equipped with ICCD camera. Maximal fluorescence signal obtained in this experiment was 23,000 RLU (Relative Light Units)..

Table 1. Comparison of EC₅₀ values of IP1-expressing Chem-1 cells.

| LIGAND | ASSAY | POTENCY (nM) | REFERENCE |
|------------------|--------------|--------------|------------------------|
| Prostaglandin I2 | Calcium Flux | 3200 | Eurofins Internal Data |

ASSAY SETUP

Fluorescence

Table 3. 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 | 95 µl (50 µl for 384-well) |
| Dispense Speed | 50 µl/sec |
| Expel Volume | 0 µl |
| Analysis | Subtract Bias Sample 1 |

Table 4. Assay Materials (Not provided)

| Description | Supplier and Product Number |
|----------------|-----------------------------|
| HBSS | Invitrogen: 14025 |
| HEPES 1M Stock | EMD Millipore: TMS-003-C |

| | |
|--|---------------------|
| Probenicid | Sigma: P8761 |
| Quest Fluo-8™, AM | AAT Bioquest: 21080 |
| Prostaglandin I2 | Various |
| Non-Binding 96/384 well Plates (for ligand prep) | Corning: 3605/ 3574 |
| Black (clear Bottom) cell assay plates | Corning: 3904/ 3712 |

HOST CELL

Chem-1, an adherent cell line expressing the promiscuous G-protein, Gα15.

EXONGENOUS GENE EXPRESSION

Human IP1 cDNA (Accession Number: NM_000960; see CODING SEQUENCE below) expressed from a proprietary E5 promoter plasmid

CODING SEQUENCE

```

atggcgggattcgtgcaggaacctcacctacgtgcggggctcgggtggggcgggccaccagc
M A D S C R N L T Y V R G S V G P A T S
accctgatgttcgtggccggtgtggtgggcaacgggctggccctgggcatcctgagcgca
T L M F V A G V V G N G L A L G I L S A
cggcgaccggcgcgccctcggccttcgcggtgctggtcaccggactggcggccaccgac
R R P A R P S A F A V L V T G L A A T D
ctgctgggaccagcttcctgagcccgccgtgttcgtggcctatgcgcgcaacagctcc
L L G T S F L S P A V F V A Y A R N S S
ctgctgggctggcccgaggcgcccccgcctgtgcatgcttcgcttcgcatgacc
L L G L A R G G P A L C D A F A F A M T
ttcttcggcctggcgtccatgctcctccttggcctggcctgggagcgtgctgctggcg
F F G L A S M L I L F A M A V E R C L A
ctgagccaccctacctctacgcgcagctggacgggccccgctgccccgctggcgctg
L S H P Y L Y A Q L D G P R C A R L A L
ccagccatctacgccttctgctccttctgcgcgctgccccctgctgggctgggccc
P A I Y A F C V L F C A L P L L G L G Q
caccagcagtaactgccccggcagctggtgcttcctcccgatgctgctgggcccagccggg
H Q Q Y C P G S W C F L R M R W A Q P G
ggcggcgccttctcgtggcctacgcccggcctgggtggccctgctggtggctgccccttc
G A A F S L A Y A G L V A L L V A A I F
ctctgcaacggctcggtcaccctcagcctctgcccgatgtaccgccagcagaagcgccac
L C N G S V T L S L C R M Y R Q Q K R H
cagggctcctcgggtccacggcgcgaccggagaggacgaggtggaccacctgatcctg
Q G S L G P R P R T G E D E V D H L I L
ctggccctcatgacagtggctcatggcgtgtgctccctgcctctcagatccgctgcttc
L A L M T V M A V C S L P L T I R C F
accaggtgctgccccctgacagcagcagtgagatgggggacctccttgccttccgcttc
T Q A V A P D S S S E M G D L L A F R F
tacgccttcaacccatcctggaccctgggtcttccatcctttccgcaaggctgtcttc
Y A F N P I L D P W V F I L F R K A V F
cagcgactcaagctcctgggtctgctgctgctgctcgggctgcccacggagactcgcag
Q R L K L W V C C L C L G P A H G D S Q
acacccctttccagctcgcctccgggaggaggacccaaggccccctctgctcctgtg
T P L S Q L A S G R R D P R A P S A P V
ggaaaggaggggagctgctgcttcttgcgcttggggcgaggggcagggtggagcccttg
G K E G S C V P L S A W G E G Q V E P L
cctcccacacagcagctccagcggcagcgcctgggaacgtcgtccaagcagaagccagc
P P T Q Q S S G S A V G T S S K A E A S
gtcgcctgctccctctgctga
V A C S L C -

```

RELATED PRODUCTS

PRODUCT NUMBER

DESCRIPTION

HTSCHEM-1

ChemiScreen™ Chem-1 Parental Cell Line (control cells)

REFERENCES

1. Narumiya S, Sugimoto Y and Ushikubi F (1999) Prostanoid receptors: structures, properties, and functions. *Physiol. Rev.* 79: 1193–1226.
2. Murata T, Ushikubi F, Matsuoka T et al. (1997) Altered pain perception and inflammatory response in mice lacking prostacyclin receptor, *Nature* 388: 678–682.
3. Cheng Y, Austin SC, Rocca B et al. (2002) Role of prostacyclin in the cardiovascular response to thromboxane A2. *Science* 296: 539–541.
4. Xiao CH, Hara A, Yuhki KI et al. (2001) Roles of prostaglandin I2 and thromboxane A2 in cardiac ischemia-reperfusion injury: a study using mice lacking their respective receptors, *Circulation* 104: 2210–2215.
5. Fitzgerald GA (2004) Coxibs and cardiovascular disease. *N. Engl. J. Med.* 351: 1709–1711.

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 **this product**

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