

PRODUCT DATASHEET

**Ready-to-Assay™ ChemiBrite™ IP1
Prostanoid Receptor Frozen Cells****CATALOG NUMBER: HTS131LRTA****CONTENTS:** Pack contains 2 vials of mycoplasma-free cells, 1 ml per vial.**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 overnight recovery, assays for calcium response.

ChemiBrite™ cells express a novel variant of clytin, a calcium-activated photoprotein, to enable sensitive luminescent detection of ligand-induced calcium flux. The ChemiBrite™ version of clytin contains a mutation that increases its affinity for calcium to a level that permits detection of cytosolic calcium in many cells with greater sensitivity than other photoproteins targeted to the mitochondria. Luminescent calcium assays offer several advantages over fluorescent calcium assays including increased sensitivity and lack of interference from fluorescent compounds.

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). Cloned human IP₁ receptor-expressing ChemiBrite™ cells are made by stable transfection of HEK293 cells with ChemiBrite™ clytin and IP₁ Receptor. These stability tested cells are ready for luminescent analysis of agonists, antagonists and modulators at the IP₁ receptor.

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: Luminescent Mode and Fluorescent Mode ; cAMP accumulation

APPLICATION DATA

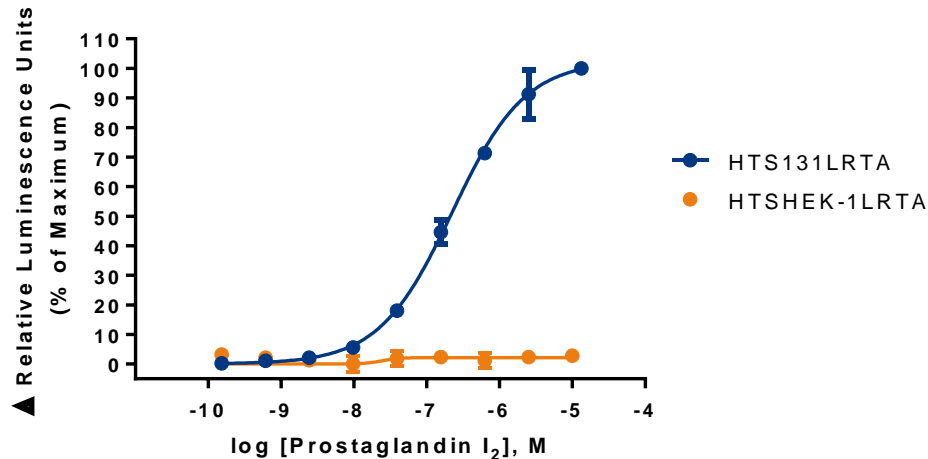


Figure 1. Representative data for activation of IP1 receptor stably expressed in HEK293 cells induced by Prostaglandin I₂ using a luminescent calcium flux assay. IP1-expressing HEK293 cells were loaded with 10 μM coelenterazine for 3 hrs at room temperature and calcium flux in response to the indicated ligand(s) was determined on a Molecular Devices FLIPRTETRA® with ICCD camera in 96-well format with a final concentration of 0.5% DMSO. Luminescence signal obtained in this experiment was 30,000 RLU (Relative Light Units) as measured by AUC (area under curve) for 80 sec post agonist addition using the provided protocol. Similarly parental cells (catalog #: HTSHEK-1L) were tested to determine the specificity of the resulting signal.

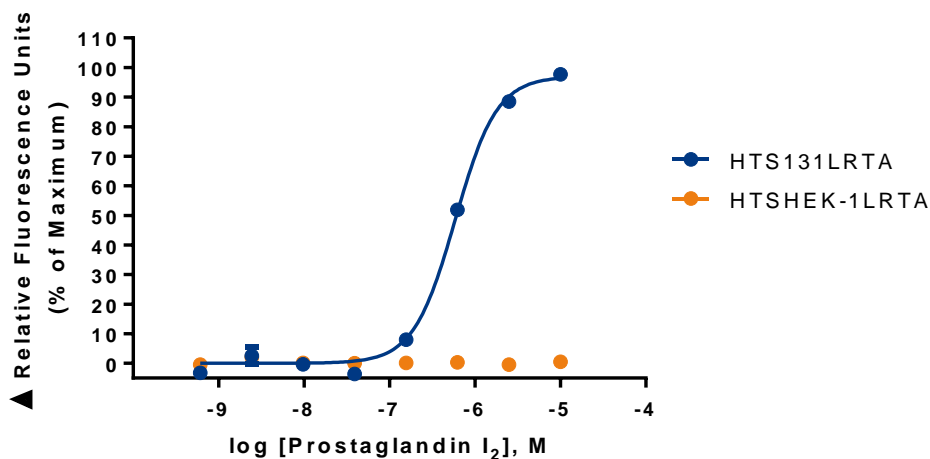


Figure 2. Representative data for activation of IP1 receptor stably expressed in HEK293 cells induced by Prostaglandin I₂ using a fluorescent calcium flux assay. IP1-expressing HEK293 cells were seeded at 50,000 cells per well into a 96-well plate. The following day the cells were loaded with a calcium dye, and calcium flux in response to the indicated ligand(s) with a final concentration of 0.5% DMSO was determined on a Molecular Devices FLIPRTETRA® with ICCD camera. Maximal fluorescence signal obtained in this experiment was 3,000 RLU. Similarly parental cells (catalog #: HTSHEK-1L) were tested to determine the specificity of the resulting signal.

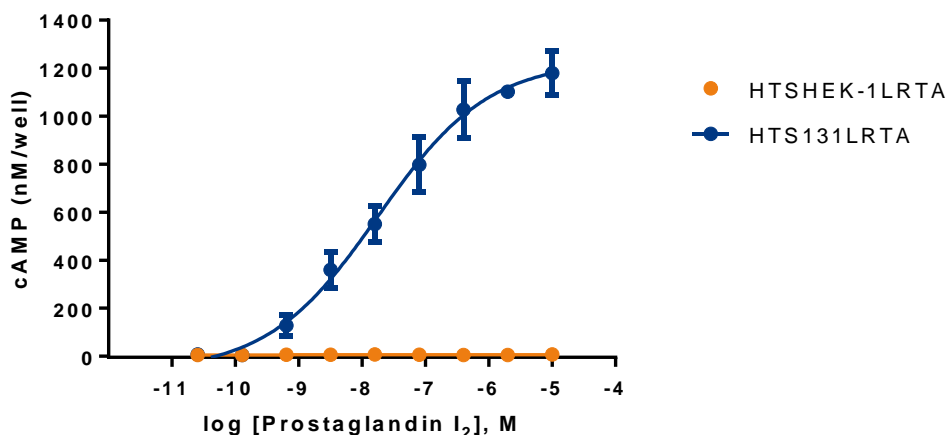


Figure 3. Representative data for activation of IP₁ receptor stably expressed in HEK293 cells induced by Prostaglandin I₂ using a cAMP accumulation assay. IP₁-expressing HEK293 cells were seeded at 25,000 cells per well into a 96-well plate, and the following day the cells were treated with Prostaglandin I₂ for 10 minutes in the presence of 100 μ M IBMX and 0.5% DMSO to determine receptor mediated cAMP generation using a time-resolve fluorescence resonance energy transfer (TR-FRET) assay measured on the BioTek Synergy. Similarly parental cells (catalog #: HTSHEK-1L) were tested to determine the specificity of the resulting signal.

Table 1. EC₅₀ values of IP₁-expressing HEK293 cells

LIGAND	ASSAY	POTENCY (nM)	REFERENCE
Prostaglandin I ₂	Calcium Flux - Luminescence	222	Eurofins Internal Data
Prostaglandin I ₂	Calcium Flux - Fluorescence	590	Eurofins Internal Data
Prostaglandin I ₂	cAMP accumulation	18	Eurofins Internal Data

ASSAY SETUP

Luminescence

Table 2. Settings for FLIPR^{TETRA}® with ICCD camera option

Option	Setting
Read Mode	Luminescence
Ex/Em	None/None
Camera Gain	280,000
Gate Open	100 %
Exposure Time	0.9 sec
Read Interval	1 sec.
Dispense Volume	50 μ l (25 μ l for 384-well)
Dispense Height	95 μ l (50 μ l for 384-well)
Dispense Speed	50 μ l/sec
Expel Volume	0 μ l
Analysis	Subtract Bias Sample 1

Table 3. Luminescence Assay Materials (Not provided)

Description	Supplier and Product Number
HBSS	Invitrogen: 14025
HEPES 1M Stock	Millipore Sigma: H3537
Quest Fluo-8 TM , AM	AAT Bioquest: 21080
Prostaglandin I ₂ ligand	Cayman: 18220

Non-Binding 96/384 well Plates (for ligand prep)	Corning: 3605/ 3574
Black (clear Bottom) cell assay plates	Corning: 3904/ 3712
Coelenterazine-h (250µg). Prepare to 10mM	Promega: S2011

Fluorescence

Table 4. 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 5. Fluorescence Assay Materials (Not provided)

Description	Supplier and Product Number
HBSS	Invitrogen: 14025
HEPES 1M Stock	Millipore Sigma: H3537
Probenicid	Sigma: P8761
Quest Fluo-8 TM , AM	AAT Bioquest: 21080
Prostaglandin I2 ligand	Cayman: 18220
Non-Binding 96/384 well Plates (for ligand prep)	Corning: 3605/ 3574
Black (clear Bottom) cell assay plates	Corning: 3904/ 3712

cAMP

Table 6. Settings for Plate Reader

Option	Setting
Excitation	300 nm
Emission	665/620 nm

Table 7. cAMP Assay Materials (Not provided)

Description	Supplier and Product Number
HEPES 1M Stock	Millipore Sigma: H3537
IBMX Buffer	Sigma #I5879
96-Well Flat Bottom Microtiter Plates	Costar #3917
Non-Binding 96 well Plates (for ligand prep)	Costar: #3789
Prostaglandin I2 ligand	Cayman: 18220
cAMP Hi Range Kit	CisBio # 62AM6PEC

Assay Protocol – Luminescence

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 18-24 h.
9. Next day, prepare Assay buffer (HBSS, 20mM HEPES, pH 7.4) and Loading buffer (Assay buffer with 10 μ M coelenterazine). *Note: Please prepare coelenterazine stock according to Manufacturer's Recommendations at 10mM to allow for 1:1000 dilution into Loading buffer (10 μ M final concentration). It is critical that coelenterazine solution is prepared at room temperature and is protected from light.*
10. Remove plate from incubator and quickly invert plate on an absorbent pad and blot to remove all Media Component.
11. Add Loading buffer to assay plate (100 μ L/well for 96-well plate). Incubate plate for 1.5 h at room temperature, protected from light.
12. Prepare ligands in assay buffer at 3x final concentration in non-binding plates. Use Buffer Only Control Wells for Background Subtraction.
13. Create protocol for ligand addition. Please refer to FLIPR^{TETRA}® settings provided in Table 2. Set time course for 180 s, with ligand addition at 10 s.
14. After the run is complete, apply subtract bias on sample 1. Export data to analyze using the area under the curve statistic.

Assay Protocol – Fluorescence

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. Move assay plate to a humidified 37°C 5% CO₂ incubator for 18-24 h.
8. Next day, prepare Assay buffer (HBSS, 20mM HEPES, 2.5 mM Probenicid, pH 7.4) and Loading buffer (Assay buffer with 5 mM Fluo8 Dye). *Note: Please prepare Fluo8 stock according to Manufacturer's Recommendations*
9. Remove plate from incubator and quickly invert plate on an absorbent pad and blot to remove all Media Component.
10. Add Loading buffer to assay plate (100 μ L/well for 96-well plate). Incubate plate for 1.5 h at room temperature, protected from light.
11. Prepare ligands in assay buffer at 3x final concentration in non-binding plates. Use Buffer Only Control Wells for Background Subtraction.
12. Create protocol for ligand addition. Please refer to FLIPR^{TETRA}® settings provided in Table 2. Set time course for 180 s, with ligand addition at 10 s.
13. After the run is complete, apply subtract bias on sample 1. We recommend using Negative Control Correction with Buffer Only Wells. Export data to according to research needs. For most Calcium Flux analysis using Export of Max Signal to end of run is sufficient.

Assay Protocol – cAMP

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 20 mL of pre-warmed Media Component to resuspend the cell pellet.
6. Seed cell suspension into appropriate assay microplate (100 µL/well for 96-well plate, 25 µL/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 18-24 h.
9. Prepare Assay Buffer (HBSS, 20mM HEPES, pH 7.4)
10. Prepare anti-cAMP-Cryptate and cAMP-D2 conjugate stocks. Please prepare stocks according to the Manufacturers Recommendations for reconstitution volume to be used according to Appendix-2 of the insert (Two-step protocol). Mix gently. Store at 10 to 35 °C after reconstitution. Reagent can only go through one cycle of freeze and thaw.
11. Prepare 1M IBMX Buffer: add 450 µL of DMSO to 1 vial of 100 mg IBMX powder.
12. Next day, prepare 2mM IBMX Solution: Place the 1M IBMX Buffer and the Assay Buffer in 50-60°C water bath for 10 min to 15 min. Add 10 µL of 1M IBMX Buffer to 4.990 mL of Assay Buffer. Mix by vortex. Place the 2 mM of IBMX Buffer into 34-40°C Incubator for at least 10 minutes or until ready to add to cells.
13. Prepare 25uL/well of compounds for testing.
14. Remove the cell assay plate from previous day from cell culture incubator. Quickly invert plate on an absorbent pad and blot to remove all Media Component. Add 25 µL of the 2 mM IBMX Buffer to the plate with the seeded GLP-1/HEK cells. Tap plate gently 3-4 times. Cover plate and incubate inside 34-40°C incubator, static until ready for compound addition.
15. Add 25 µL compounds, internal control and test sample dilutions to cell assay plate. Cover plate and incubate for 15 to 25 min at 20-25°C
16. Prepare fresh working dilutions of 1:24 of anti-cAMP-Cryptate and cAMP-D2 conjugate in Lysis Buffer. Protect from light. Do not vortex.
17. Add 125 µL of cAMP-D2 solution into 2,875 µL of Lysis Buffer for total volume of 3 mL.
18. Add 125 µL of anti-cAMP-Cryptate solution into 2,875 µL of Lysis Buffer for total volume of 3 mL.
19. It is imperative that detection reagents are added to plate in the following order:
Add 25 µL cAMP-D2 conjugate/lysis buffer solution to each well of cell assay plate.
Add 25 µL anti-cAMP-Cryptate/lysis buffer solution to each well of cell assay plate.
20. Cover with aluminum plate sealer and incubate the cell assay plate 20-25°C, for 60 to 65 min (If available, use gentle plate shaker). Protect from light.
21. Read plate on a plate reader with 330 nm (excitation) and 665/620 nm (emission).
22. Calculate Ratio Emission 665/620 nm.

HOST CELL

HEK293

EXOGENOUS GENE EXPRESSION

PTGIR cDNA (Accession Number: NM_000960; see CODING SEQUENCE below) and a proprietary mutant clytin photoprotein each expressed in a bicistronic vector

CODING SEQUENCE

```

ATG GCG GAT TCG TGC AGG AAC CTC ACC TAC GTG CGG GGC TCG GTG GGG CCG GCC
M  A  D  S  C  R  N  L  T  Y  V  R  G  S  V  G  P  A

ACC AGC ACC CTG ATG TTC GTG GCC GGT GTG GTG GGC AAC GGG CTG GCC CTG GCC
T  S  T  L  M  F  V  A  G  V  V  G  N  G  L  A  L  G

```

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

RELATED PRODUCTS

PRODUCT NUMBER	DESCRIPTION
HTSHEK-1L	ChemiBrite™ HEK stable cell line (control cells)
HTS081L	ChemiBrite™ TP Prostanoid family Stable Cell Line
HTS081LRTA	Ready-to-Assay™ ChemiBrite™ TP Prostanoid Receptor Frozen Cells
HTS185L	ChemiBrite™ HEK stable EP ₂ Prostanoid Receptor Cell Line
HTS185LRTA	Ready-to-Assay™ ChemiBrite™ EP ₂ Prostanoid Receptor Frozen Cells
HTS142L	ChemiBrite™ HEK stable EP ₄ Prostanoid Receptor Cell Line
HTS142LRTA	Ready-to-Assay™ ChemiBrite™ EP ₄ Prostanoid Receptor Frozen Cells
HTS091L	ChemiBrite™ HEK stable DP Prostanoid Receptor Cell Line
HTS091LRTA	Ready-to-Assay™ ChemiBrite™ DP Prostanoid Receptor Frozen Cells
HTS092L	ChemiBrite™ HEK stable EP ₃ Prostanoid Receptor Cell Line
HTS092LRTA	Ready-to-Assay™ ChemiBrite™ EP ₃ Prostanoid Receptor Frozen Cells
HTS093L	ChemiBrite™ HEK stable FP Prostanoid Receptor Cell Line
HTS093LRTA	Ready-to-Assay™ ChemiBrite™ FP Prostanoid Receptor Frozen Cells
HTS099L	ChemiBrite™ HEK stable EP ₁ Prostanoid Receptor Cell Line
HTS099LRTA	Ready-to-Assay™ ChemiBrite™ EP ₁ Prostanoid Receptor Frozen Cells
HTS131M	ChemiScreen™ IP1 Prostanoid receptor membrane prep
HTS131RTA	Ready-to-Assay™ IP1 Prostanoid receptor frozen cells
HTSCHEM-1RTA	Ready-to-Assay™ Chem-1 host frozen cells (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 **Ready-to-Assay™ ChemiBrite™ IP1 Prostanoid Receptor Frozen Cells, Product No. HTS131LRTA**

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.