

PRODUCT DATASHEET**Ready-to-Assay™ GPBA Bile Acid
Receptor Frozen Cells****CATALOG NUMBER: HTS238LRTA**

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.

ChemiBrite cells co-express a GPCR along with 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 mitochondrially expressed photoproteins. Luminescent calcium assays offer several advantages over fluorescent calcium assays including; lower substrate cost, increased sensitivity, and lack of interference from fluorescent compounds.

The bile acid lithocholic acid acts as an agonist on the GPBA GPCR, also known as GPR131 or TGR5, to increase intracellular cAMP in cells that express the receptor (Maruyama et al., 2002). Bile acids and other GPBA ligands increase glucagon-like peptide-1 (GLP-1) secretion from intestinal cells. Although Type II diabetics have impaired ability to secrete GLP-1, their responsiveness to this compound, which enhances the effects of insulin, is normal. Thus GPBA agonists are of interest in developing therapeutics for diabetes. Interestingly, it has been suggested that the anti-diabetic effect of olive leaves could in part be mediated by agonist activity at GPBA receptors by oleanolic acid (Sato et al. 2007). Cloned GPBA receptor-expressing ChemiBrite cells were made by stable transfection of HEK293 cells with ChemiBrite clytin and the receptor and a promiscuous G protein to couple the receptor to the calcium signaling pathway. The cells have been cryopreserved at an optimal time post-transfection. Upon thaw, recovery, and loading, the cells are ready for luminescent, fluorescent and cAMP accumulation analysis of agonists, antagonists and modulators at the GPBA 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.

**Eurofins Pharma Bioanalytics
Services US Inc.**

15 Research Park Drive
St Charles MO 63304
USA

T | +1 844 522 7787
F | +1 636 362 7131
www.eurofins.com

APPLICATIONS

Calcium Flux Assays: Luminescent Mode and Fluorescent Mode; cAMP accumulation

APPLICATION DATA

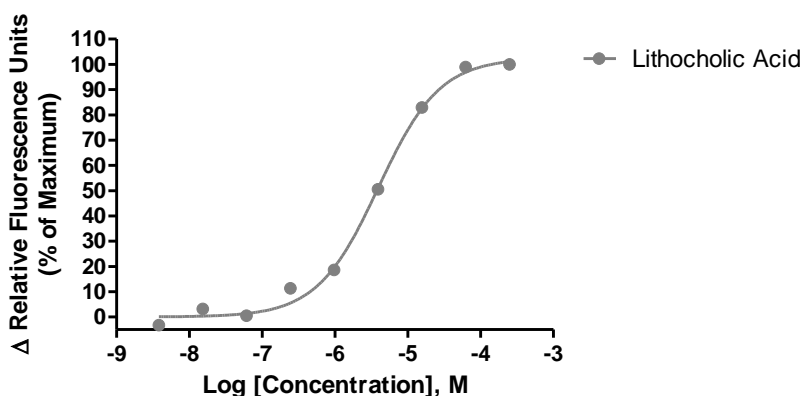


Figure 1. Representative data for activation of GPBA receptor. Calcium flux in GPBA-expressing HEK293 cell line induced by Lithocholic Acid. GPBA-expressing HEK293 cells were loaded with a calcium dye, and calcium flux in response to the indicated ligand(s), 4-fold serial dilution with each concentration performed in triplicate, was determined on a Molecular Devices FLIPR^{TETRA}. Maximal fluorescence signal obtained in this experiment was 8,000 RLU (Relative Light Units).

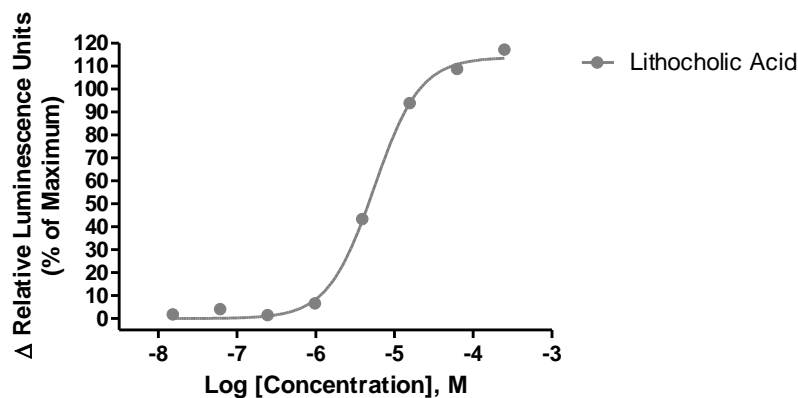


Figure 2. Representative data for activation of GPBA receptor expressed in HEK293 cells induced by Lithocholic Acid using a luminescent calcium flux assay. GPBA-expressing HEK293 cells were loaded with 10 μ M coelenterazine for 2h at room temperature and calcium flux in response to the indicated ligand was determined on a Molecular Devices FLIPR^{TETRA}® with ICCD camera in 96-well format. Luminescence signal obtained in this experiment was 35,000 RLU (Relative Light Units) as measured by area-under-curve for 80s post agonist addition using the provided protocol.

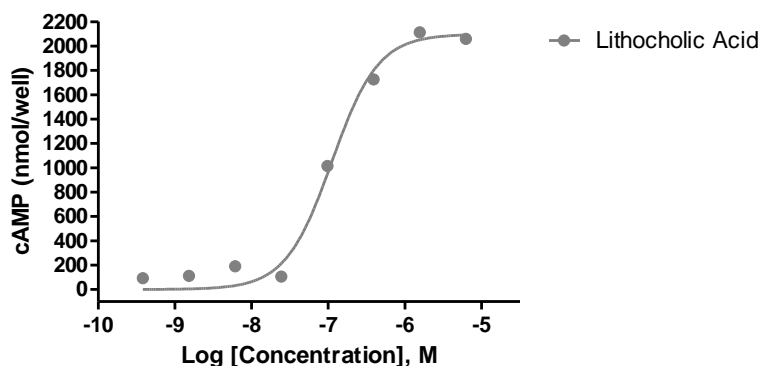


Figure 3. Representative data for activation of GPBA receptor stably expressed in HEK293 cells induced by Lithocholic Acid using a cAMP accumulation assay. GPBA-expressing HEK293 cells were seeded into a 96-well plate, and the following day the cells were treated with Lithocholic Acid for 15 minutes in the presence of 100 μ M IBMX to determine receptor-mediated cAMP generation using a time-resolved fluorescence resonance energy transfer (TR-FRET) assay measured on the BioTek Synergy.

Table 1. Comparison of EC₅₀ values of GPBA-expressing HEK293 cells.

| LIGAND | ASSAY | POTENCY (μ M) | REFERENCE |
|------------------|-----------------------------|--------------------|------------------------|
| Lithocholic Acid | Calcium Flux - Fluorescence | 3.9 | Eurofins Internal Data |
| Lithocholic Acid | Calcium Flux - Luminescence | 5.5 | Eurofins Internal Data |
| Lithocholic Acid | cAMP accumulation | 0.11 | 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 |

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 TM , AM | AAT Bioquest: 21080 |
| Lithocholic Acid ligand | Sigma: T7515 |
| 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 |

Assay Protocol – Luminescence

1. Dissociate Culture as Recommended. Collect in Basal Medium. Document Cell Count and Viability
2. Centrifuge the cell suspension at 190 x g for six min
3. Remove supernatant. Gently resuspend the cell pellet in Basal Medium. *It is suggested that end user optimize cell plating based on individual formats.* (Default: Resuspend in volume to achieve 5×10^5 cells/ml (i.e, if collected 5×10^6 TC, $\frac{5 \times 10^6}{5 \times 10^5/ml} = 10$ mL volume)
4. Seed cell suspension into black, clear bottom plate (100 µL/well for 96-well plate). *When seeding is complete, place the assay plate at room temperature for 30 min.*
5. Move assay plate to a humidified 37°C 5% CO₂ incubator for 18-24 h.
6. Next day, prepare Assay buffer (HBSS, 20mM HEPES, pH 7.4) and Loading buffer (Assay buffer with 10µM coelenterazine). *Note: Please prepare coelenterazine stock according to Manufacturer's Recommendations at 10mM to allow for 1:1000 dilution into Loading buffer (10uM final concentration). It is critical that coelenterazine solution is prepared at room temperature and is protected from light.*
7. Remove medium from assay plate and wash 1X with Assay Buffer.
8. Add Loading buffer to assay plate (100 µL/well for 96-well plate). Incubate plate for 1.5 h at room temperature, protected from light.
9. Prepare ligands in assay buffer at 3x final concentration in non-binding plates. Use Buffer Only Control Wells for Background Subtraction.
10. 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.
11. 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. Dissociate Culture as Recommended. Collect in Basal Medium. Document Cell Count and Viability
2. Centrifuge the cell suspension at 190 x g for six min
3. Remove supernatant. Gently resuspend the cell pellet in Basal Medium. *It is suggested that end user optimize cell plating based on individual formats.* (Default: Resuspend in volume to achieve 5×10^5 cells/ml (i.e, if collected 5×10^6 TC, $\frac{5 \times 10^6}{5 \times 10^5/ml} = 10$ mL volume)
4. Seed cell suspension into black, clear bottom plate (100 μ L/well for 96-well plate). *When seeding is complete, place the assay plate at room temperature for 30 min.*
5. Move assay plate to a humidified 37°C 5% CO₂ incubator for 18-24 h.
6. 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*
7. Remove medium from assay plate and wash 1X with Assay Buffer.
8. 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.
9. Prepare ligands in assay buffer at 3x final concentration in non-binding plates. Use Buffer Only Control Wells for Background Subtraction.
10. 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.
11. 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.

HOST CELL

HEK293

EXOGENOUS GENE EXPRESSION

GPBAR1 cDNA (Accession Number: NM_170699; see CODING SEQUENCE below) and a proprietary mutant clytin photoprotein, and promiscuous G protein each expressed in a bicistronic vector

CODING SEQUENCE

```

ATG ACG CCC AAC AGC ACT GGC GAG GTG CCC AGC CCC ATT CCC AAG GGG GCT TTG
M T P N S T G E V P S P I P K G A L

GGG CTC TCC CTG GCC CTG GCA AGC CTC ATC ATC ACC GCG AAC CTG CTC CTA GCC
G L S L A L A S L I I T A N L L L A

CTG GGC ATC GCC TGG GAC CGC CGC CTG CGC AGC CCA CCT GCT GGC TGC TTC TTC
L G I A W D R R L R S P P A G C F F

CTG AGC CTA CTG CTG GCT GGG CTG CTC ACG GGT CTG GCA TTG CCC ACA TTG CCA
L S L L L A G L L T G L A L P T L P

GGG CTG TGG AAC CAG AGT CGC CGG GGT TAC TGG TCC TGC CTC CTC GTC TAC TTG
G L W N Q S R R G Y W S C L L V Y L

GCT CCC AAC TTC TCC TTC CTC TCC CTG CTT GCC AAC CTC TTG CTG GTG CAC GGG
A P N F S F L S L L A N L L L V H G

GAG CGC TAC ATG GCA GTC CTG AGG CCA CTC CAG CCC CCT GGG AGC ATT CGG CTG
E R Y M A V L R P L Q P P G S I R L

GCC CTG CTC CTC ACC TGG GCT GGT CCC CTG CTC TTT GCC AGT CTG CCC GCT CTG
A L L L T W A G P L L F A S L P A L

GGG TGG AAC CAC TGG ACC CCT GGT GCC AAC TGC AGC TCC CAG GCT ATC TTC CCA
G W N H W T P G A N C S S Q A I F P

GCC CCC TAC CTG TAC CTC GAA GTC TAT GGG CTC CTG CTG CCC GCC GTG GGT GCT
A P Y L Y L E V Y G L L L P A V G A
    
```

```

GCT GCC TTC CTC TCT GTC CGC GTG CTG GCC ACT GCC CAC CGC CAG CTG CAG GAC
A A F L S V R V L A T A H R Q L Q D

ATC TGC CGG CTG GAG CGG GCA GTG TGC CGC GAT GAG CCC TCC GCC CTG GCC CGG
I C R L E R A V C R D E P S A L A R

GCC CTT ACC TGG AGG CAG GCA AGG GCA CAG GCT GGA GCC ATG CTG CTC TTC GGG
A L T W R Q A R A Q A G A M L L F G

CTG TGC TGG GGG CCC TAC GTG GCC ACA CTG CTC CTC TCA GTC CTG GCC TAT GAG
L C W G P Y V A T L L L S V L A Y E

CAG CGC CCG CCA CTG GGG CCT GGG ACA CTG TTG TCC CTC CTC TCC CTA GGA AGT
Q R P P L G P G T L L S L L S L G S

GCC AGT GCA GCG GCA GTG CCC GTA GCC ATG GGG CTG GGC GAT CAG CGC TAC ACA
A S A A A V P V A M G L G D Q R Y T

GCC CCC TGG AGG GCA GCC GCC CAA AGG TGC CTG CAG GGG CTG TGG GGA AGA GCC
A P W R A A A Q R C L Q G L W G R A

TCC CGG GAC AGT CCC GGC CCC AGC ATT GCC TAC CAC CCA AGC AGC CAA AGC AGT
S R D S P G P S I A Y H P S S Q S S

GTC GAC CTG GAC TTG AAC TAA
V D L D L N STP

```

RELATED PRODUCTS

PRODUCT NUMBER

HTSHEK-6L

DESCRIPTION

ChemiBrite™ HEK Parental cell line with Gα16 (control cells)

REFERENCES

1. Maruyama et al. (2002) Identification of membrane-type receptor for bile acids (M-BAR). *Biochem Biophys Res Commun* 298:714-719
2. Sato et al. (2007) Anti-hyperglycemic activity of a TGR5 agonist isolated from *Olea europaea*. *Biochem Biophys Res Commun* 362:793-798

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™ GPBA Bile Acid Receptor Frozen Cells, Product No. HTS238LRTA**

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.