

PRODUCT DATASHEET

Ready-to-Assay™ GPR109A Nicotinic Acid Receptor Frozen Cells

CATALOG NUMBER: HTS201RTA

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 overnight recovery, assays for calcium response.

Nicotinic acid (niacin), a vitamin of the B complex, is used clinically in high doses to decrease total plasma levels of cholesterol. Interestingly, the total cholesterol levels and low-density lipoprotein (LDL) concentrations decrease, while the high-density lipoprotein (HDL) concentrations increase with nicotinic acid treatment. The cholesterol-lowering effects of nicotinic acid result from inhibition of lipolysis in adipose tissue, which decreases plasma levels of free fatty acid (FFA) (Altschul et al., 1955; Carlson, 1963). In a study of nicotinic acid and myocardial infarction in the Coronary Drug Project (1966-1975), patients receiving 3 g/day nicotinic acid exhibited reduced rates of myocardial infarction (Coronary Drug Project Research Group, 1975). However, an unwanted effect of high doses of nicotinic acid is vasodilatation, occurring mainly in the upper body and face, known as flushing. Recently two receptors for nicotinic acid have been identified as class A G protein-coupled receptors that couple to Gi to inhibit accumulation of cAMP (Offermans, 2006). GPR109A (also known as HM74A in humans and PUMA-G in mice) is a high affinity receptor for nicotinic acid, whereas GPR109B (also known as HM74) is a low affinity receptor for nicotinic acid that is found in humans but not rodents (Wise et al., 2003). GPCR109A is found primarily in adipose tissue and immune cells. Cloned human GPR109A-expressing cell line is made in the Chem-4 host, which supports high levels of recombinant GPR109A expression on the cell surface and contains optimized levels of a recombinant promiscuous G protein to couple the receptor to the calcium signaling pathway. Thus, the cell line is an ideal tool for screening for agonists, antagonists, and modulators at GPR109A.

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

APPLICATION DATA

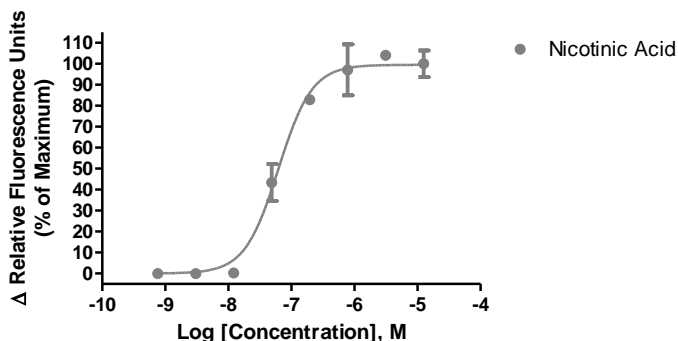


Figure 1. Representative data for activation of GPR109A receptor. Calcium flux in GPR109A –expressing Chem-4 cell line induced by Nicotinic Acid. GPR109A –expressing Chem-4 cells were loaded with a calcium dye, and calcium flux in response to the indicated ligand(s), 4-fold serial dilution with each concentration performed in duplicate, was determined on a Molecular Devices FLIPR^{TETRA} with ICCD camera. Maximal fluorescence signal obtained in this experiment was 7,500 RLU (Relative Light Units).

Table 1. EC₅₀ values of GPR109A -expressing Chem-4 cells.

LIGAND	ASSAY	POTENCY (nM)	REFERENCE
Nicotinic Acid	Calcium Flux	60	Eurofins Internal Data
Acifran (Data not shown)	Calcium Flux	250	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 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 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) Ca²⁺ dye by dissolving 1mg of Fluo-8 NW in 200 µL of DMSO. Once dissolved place 10 µL of Fluo-8 NW Ca²⁺ dye solution into 10 mL of HBSS 20mM HEPES, 2.5mM Probenecid pH 7.4 buffer and apply to assay microplate (Ca²⁺ 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^{TETRA}) or 485 nm (FLIPR1, FLIPR2, FLIPR3) and emission wavelength at 515-565 nm (FLIPR^{TETRA}) or emission filter for Ca²⁺ 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
Nicotinic Acid ligand	Sigma: N4126
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^{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	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-4, an adherent rat hematopoietic cell line expressing endogenous Gα15 protein as well as an exogenous proprietary promiscuous Gα protein.

EXONGENOUS GENE EXPRESSION

GPR109A cDNA (Accession Number: NM_177551.3; see CODING SEQUENCE below) expressed from a proprietary pHS plasmid.

CODING SEQUENCE

```

1 - ATG AAT CGG CAC CAT CTG CAG GAT CAC TTT CTG GAA ATA GAC AAG AAG AAC TGC TGT GTG TTC CGA GAT - 69
1 - M N R H H L Q D H F L E I D K K N C C V F R D - 23

70 - GAC TTC ATT GTC AAG GTG TTG CCG CCG GTG TTG GGG CTG GAG TTT ATC TTC GGG CTT CTG GGC AAT GGC - 138
24 - D F I V K V L P P V L G L E F I F G G L L G N G - 46

139 - CTT GCC CTG TGG ATT TTC TGT TTC CAC CTC AAG TCC TGG AAA TCC AGC CGG ATT TTC CTG TTC AAC CTG - 207
47 - L A L W I F C F H L K S W K S S R I F L F N L - 69

208 - GCA GTG GCT GAC TTT CTA CTG ATC ATC TGC CTG CCC TTC CTG ATG GAC AAC TAT GTG AGG CGT TGG GAC - 276
70 - A V A D F L L I I C L P F L M D N Y V R R W D - 92

277 - TGG AAG TTT GGG GAC ATC CCT TGC CGG CTG ATG CTC TTC ATG TTG GCT ATG AAC CGC CAG GGC AGC ATC - 345
93 - W K F G D I P C R L M L F M L A M N R Q G S I - 115

346 - ATC TTC CTC ACG GTG GTG GCG GTA GAC AGG TAT TTC CGG GTG GTC CAT CCC CAC CAC GCC CTG AAC AAG - 414
116 - I F L T V V A V D R Y F R V V H P H H A L N K - 138

415 - ATC TCC AAT CGG ACA GCA GCC ATC ATC TCT TGC CTT CTG TGG GGC ATC ACT ATT GGC CTG ACA GTC CAC - 483
139 - I S N R T A A I I S C L L W G I T I G L T V H - 161

484 - CTC CTG AAG AAG AAG ATG CCG ATC CAG AAT GGC GGT GCA AAT TTG TGC AGC AGC TTC AGC ATC TGC CAT - 552
162 - L L K K K M P I Q N G G A N L C S S F S I C H - 184

553 - ACC TTC CAG TGG CAC GAA GCC ATG TTC CTC CTG GAG TTC TTC CTG CCC CTG GGC ATC ATC CTG TTC TGC - 621
185 - T F Q W H E A M F L L E F F L P L G I I L F C - 207

622 - TCA GCC AGA ATT ATC TGG AGC CTG CGG CAG AGA CAA ATG GAC CGG CAT GCC AAG ATC AAG AGA GCC ATC - 690
208 - S A R I I W S L R Q R Q M D R H A K I K R A I - 230

691 - ACC TTC ATC ATG GTG GTG GCC ATC GTC TTT GTC ATC TGC TTC CTT CCC AGC GTG GTT GTG CGG ATC CGC - 759
231 - T F I M V V A I V F V I C F L P S V V V R I R - 253

760 - ATC TTC TGG CTC CTG CAC ACT TCG GGC ACG CAG AAT TGT GAA GTG TAC CGC TCG GTG GAC CTG GCG TTC - 828
254 - I F W L L H T S G T Q N C E V Y R S V D L A F - 276

829 - TTT ATC ACT CTC AGC TTC ACC TAC ATG AAC AGC ATG CTG GAC CCC GTG GTG TAC TAC TTC TCC AGC CCA - 897
277 - F I T L S F T Y M N S M L D P V V Y Y F S S P - 299

898 - TCC TTT CCC AAC TTC TTC TCC ACT TTG ATC AAC CGC TGC CTC CAG AGG AAG ATG ACA GGT GAG CCA GAT - 966
300 - S F P N F F S T L I N R C L Q R K M T G E P D - 322

967 - AAT AAC CGC AGC ACG AGC GTC GAG CTC ACA GGG GAC CCC AAC AAA ACC AGA GGC GCT CCA GAG GCG TTA - 1035
323 - N N R S T S V E L T G D P N K T R G A P E A L - 345

1036 - ATG GCC AAC TCC GGT GAG CCA TGG AGC CCC TCT TAT CTG GGC CCA ACC TCT CCT TAA
346 - M A N S G E P W S P S Y L G P T S P Stp

```

RELATED PRODUCTS

PRODUCT NUMBER

DESCRIPTION

HTSCHEM-1RTA

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

HTS201M

ChemiScreen™ GPR109A Nicotinic Acid receptor membrane prep

Note: Chem-4 cells are derived from Chem-1 cells.

REFERENCES

1. Altschul R *et al.* (1955) Influence of nicotinic acid on serum cholesterol in man. *Arch. Biochem.* 54: 558-559.
2. Carlson LA (1963) Studies on the effect of nicotinic acid on catecholamine stimulated lipolysis in adipose tissue *in vitro*. *Acta Med. Scand.* 173: 719-722.
3. Coronary Drug Project Research Group (1975) Clofibrate and niacin in coronary heart disease. *J. Am. Med. Assoc.* 231: 360-381.
4. Offermans S (2006) The nicotinic acid receptor GPR109A (HM74A or PUMA-G) as a new therapeutic target. *Trends Pharmacol. Sci.* 27: 384-390.
5. Wise A *et al.* (2003) Molecular identification of high and low affinity receptors for nicotinic acid. *J. Biol. Chem.* 278: 9869-9874.

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™ GPR109A Nicotinic Acid Receptor Frozen Cells, Product No. HTS201RTA**

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