

PRODUCT DATASHEET

Ready-to-Assay™ ChemiBrite EP₃
Prostanoid Receptor Frozen Cells

CATALOG NUMBER: HTS092LRTA

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.

Prostanoids are a series of arachidonic acid metabolites produced by the action of cyclooxygenase and subsequently by isomerases and synthases. Cells rapidly secrete prostanoids after synthesis, whereupon the prostanoids bind to a family of 8 GPCRs to exert their biological effects (Narumiya and FitzGerald, 2001). The prostaglandin PGE₂ causes pain, vasodilatation, immunosuppression of T cells, bone resorption and promotion of carcinogenesis. Four related GPCRs, EP₁, EP₂, EP₃ and EP₄, each bind to PGE₂, but the different G protein coupling status of each receptor leads to distinct biological effects. Further diversity is generated by alternative splicing; the human gene for EP₃ generates 9 alternatively-spliced mRNAs encoding 8 isoforms of EP₃ (Kotani *et al.*, 1997). These isoforms of EP₃ vary in sequence at their C-termini, and differ in their ability to couple to G_s, G_q or G_i (Kotani *et al.*, 1995). EP₃ receptor activity is required for fever induced by pyrogens, a response long attributed to prostaglandins by the antipyretic action of aspirin and other COX inhibitors (Ushikubi *et al.*, 1998). In animal models of allergy, PGE₂-mediated activation of EP₃ inhibits inflammation to counteract the allergy-promoting activity of PGD₂ (Kunikata *et al.*, 2005). Cloned EP₃ receptor-expressing ChemiBrite™ cells were made by stable transfection of HEK293 cells with ChemiBrite clytin, the EP₃ receptor and a promiscuous G protein to couple the receptor to the calcium signaling pathway. These stability tested cells are ready for luminescent analysis of agonists, antagonists and modulators at the EP₃ 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

APPLICATION DATA

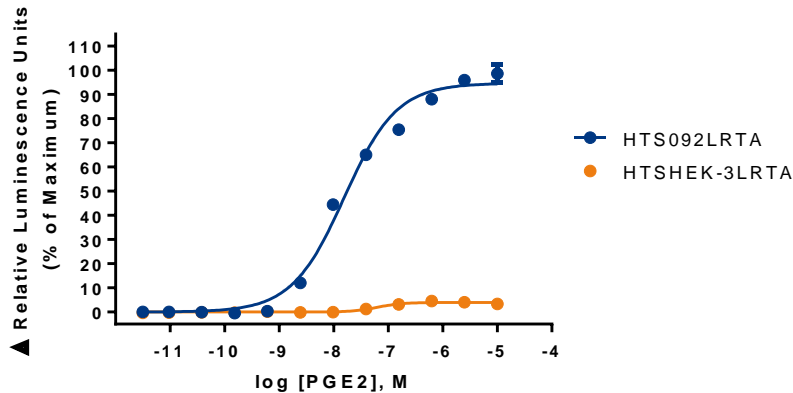


Figure 1. Representative data for activation of EP₃ receptor stably expressed in HEK293 cells induced by Prostaglandin E2 using a luminescent calcium flux assay. EP₃-expressing HEK293 cells were loaded with 10 μM coelenterazine for 3 h at room temperature and calcium flux in response to the indicated ligand was determined on a Molecular Devices FLIPR^{TETRA}® with ICCD camera in 96-well format with a final concentration of 0.5% DMSO. Luminescence signal obtained in this experiment was 200,000 RLU (Relative Light Units) as measured by AUC (area under curve) for 80 s post agonist addition using the provided protocol. Similarly parental cells (catalog #: HTSHEK-3L) were tested to determine the specificity of the resulting signal.

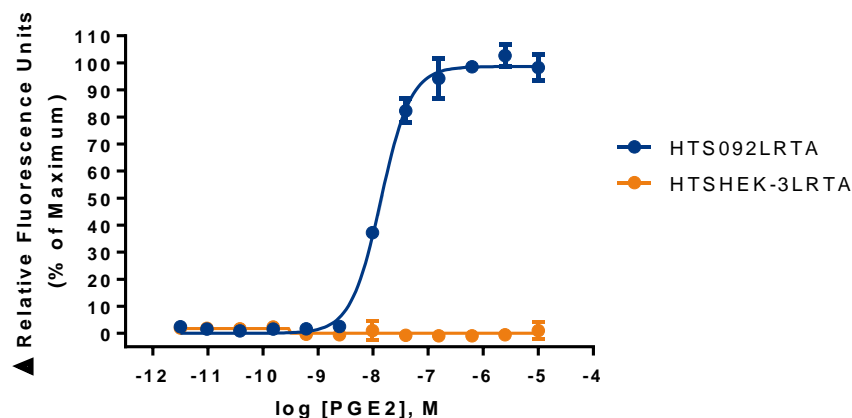


Figure 2. Representative data for activation of EP₃ receptor stably expressed in HEK293 cells induced by Prostaglandin E2 using a fluorescent calcium flux assay. EP₃-expressing HEK293 cells were seeded at 50,000 cells per well into a 96-well plate the following day the cells were loaded with a calcium dye. Calcium flux in response to the indicated ligand with a final concentration of 0.5% DMSO was determined on a Molecular Devices FLIPR^{TETRA}® with ICCD camera. Maximal fluorescence signal obtained in this experiment was 6,200 RLU as measured at 10 s post agonist addition. Similarly parental cells (catalog #: HTSHEK-3L) were tested to determine the specificity of the resulting signal.

Table 1. EC₅₀ values of EP₃-expressing HEK293 cells

| LIGAND | ASSAY | POTENCY EC ₅₀ (nM) | REFERENCE |
|------------------|-----------------------------|-------------------------------|------------------------|
| Prostaglandin E2 | Calcium Flux - Luminescence | 15 | Eurofins Internal Data |
| Prostaglandin E2 | Calcium Flux - Fluorescence | 14 | 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 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 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 |
| PGE2 ligand | Cayman: 14010 |
| 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™, AM | AAT Bioquest: 21080 |
| PGE2 ligand | Cayman: 14010 |
| Non-Binding 96/384 well Plates (for ligand prep) | Corning: 3605/ 3574 |
| Black (clear Bottom) cell assay plates | Corning: 3904/ 3712 |

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 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. 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.*
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 µL/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 µL/well for 96-well plate, 25 µL/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 µL/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.

HOST CELL

HEK293

EXONGENOUS GENE EXPRESSION

Human PTGER3 cDNA (Accession Number: NM_198716; see CODING SEQUENCE below) and a proprietary mutant clytin photoprotein and promiscuous G protein expressed in a bicistronic vector

CODING SEQUENCE

```

ATG AAG GAG ACC CGG GGC TAC GGA GGG GAT GCC CCC TTC TGC ACC CGC CTC AAC CAC TCC TAC ACA GGC
M K E T R G Y G G D A P F C T R L N H S Y T G

ATG TGG GCG CCC GAG CGT TCC GCC GAG GCG CGG GGC AAC CTC ACG CGC CCT CCA GGG TCT GGC GAG GAT
M W A P E R S A E A R G N L T R P P G S G E D

TGC GGA TCG GTG TCC GTG GCC TTC CCG ATC ACC ATG CTG CTC ACT GGT TTC GTG GGC AAC GCA CTG GCC
C G S V S V A F P I T M L L T G F V G N A L A

ATG CTG CTC GTG TCG CGC AGC TAC CGG CGC CGG GAG AGC AAG CGC AAG AAG TCC TTC CTG CTG TGC ATC
M L L V S R S Y R R R E S K R K K S F L L C I

GGC TGG CTG GCG CTC ACC GAC CTG GTC GGG CAG CTT CTC ACC ACC CCG GTC GTC ATC GTC GTG TAC CTG
G W L A L T D L V G Q L L T T P V V I V V Y L

TCC AAG CAG CGT TGG GAG CAC ATC GAC CCG TCG GGG CGG CTC TGC ACC TTT TTC GGG CTG ACC ATG ACT
S K Q R W E H I D P S G R L C T F F G L T M T

GTT TTC GGG CTC TCC TCG TTG TTC ATC GCC AGC GCC ATG GCC GTC GAG CGG GCG CTG GCC ATC AGG GCG
V F G L S S L F I A S A M A V E R A L A I R A

CCG CAC TGG TAT GCG AGC CAC ATG AAG ACG CGT GCC ACC CGC GCT GTG CTG CTC GGC GTG TGG CTG GCC
P H W Y A S H M K T R A T R A V L L G V W L A

GTG CTC GCC TTC GCC CTG CTG CCG GTG CTG GGC GTG GGC CAG TAC ACC GTC CAG TGG CCC GGG ACG TGG
V L A F A L L P V L G V G Q Y T V Q W P G T W

TGC TTC ATC AGC ACC GGG CGA GGG GGC AAC GGG ACT AGC TCT TCG CAT AAC TGG GGC AAC CTT TTC TTC
C F I S T G R G G N G T S S S H N W G N L F F

GCC TCT GCC TTT GCC TTC CTG GGG CTC TTG GCG CTG ACA GTC ACC TTT TCC TGC AAC CTG GCC ACC ATT
A S A F A F L G L L A L T V T F S C N L A T I

AAG GCC CTG GTG TCC CGC TGC CGG GCC AAG GCC ACG GCA TCT CAG TCC AGT GCC CAG TGG GGC CGC ATC
K A L V S R C R A K A T A S Q S S A Q W G R I

ACG ACC GAG ACG GCC ATT CAG CTT ATG GGG ATC ATG TGC GTG CTG TCG GTC TGC TGG TCT CCG CTC CTG
T T E T A I Q L M G I M C V L S V C W S P L L

ATA ATG ATG TTG AAA ATG ATC TTC AAT CAG ACA TCA GTT GAG CAC TGC AAG ACA CAC ACG GAG AAG CAG
I M M L K M I F N Q T S V E H C K T H T E K Q

AAA GAA TGC AAC TTC TTC TTA ATA GCT GTT CGC CTG GCT TCA CTG AAC CAG ATC TTG GAT CCT TGG GTT
K E C N F F L I A V R L A S L N Q I L D P W V

TAC CTG CTG TTA AGA AAG ATC CTT CTT CGA AAG TTT TGC CAG ATG AGA AAA AGA AGA CTC AGA GAG CAA
Y L L L R K I L L R K F C Q M R K R R L R E Q

GAG GAA TTT TGG GGA AAT TGA
E E F W G N Stp

```

RELATED PRODUCTS

| Product Number | Description |
|----------------|--|
| HTSHEK-1L | ChemiBrite™ HEK stable cell line (control cells) |
| HTSHEK-3L | ChemiBrite™ HEK Gαq stable cell line (control cells) |
| HTS185L | ChemiBrite™ HEK stable EP2 Prostanoid Receptor Cell Line |
| HTS185LRTA | Ready-to-Assay™ ChemiBrite™ EP2 Prostanoid Receptor Frozen Cells |
| HTS142L | ChemiBrite™ HEK stable EP4 Prostanoid Receptor Cell Line |
| HTS142LRTA | Ready-to-Assay™ ChemiBrite™ EP4 Prostanoid Receptor Frozen Cells |
| HTS081L | ChemiBrite™ HEK stable TP Prostanoid Receptor Cell Line |
| HTS081LRTA | Ready-to-Assay™ ChemiBrite™ TP 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 EP3 Prostanoid Receptor Cell Line |
| HTS093L | ChemiBrite™ HEK stable FP Prostanoid Receptor Cell Line |
| HTS093LRTA | Ready-to-Assay™ ChemiBrite™ FP Prostanoid Receptor Frozen Cells |
| HTS092RTA | Ready-to-Assay™ EP3 Prostanoid receptor frozen cells |
| HTS092M | ChemiScreen™ EP3 Prostanoid receptor Membrane Prep |
| HTS099L | ChemiBrite™ HEK stable EP1 Prostanoid Receptor Cell Line |
| HTS099LRTA | Ready-to-Assay™ ChemiBrite™ EP ₁ Prostanoid Receptor Frozen Cells |
| HTS131L | ChemiBrite™ HEK stable IP1 Prostanoid Receptor Cell Line |
| HTS131LRTA | Ready-to-Assay™ ChemiBrite™ IP1 Prostanoid Receptor Frozen Cells |

REFERENCES

1. Kotani M *et al.* (1995) Molecular cloning and expression of multiple isoforms of human prostaglandin E receptor EP₃ subtype generated by alternative messenger RNA splicing: multiple second messenger systems and tissue-specific distributions. *Mol Pharmacol.* 48: 869-879.
2. Kotani M *et al.* (1997) Structural Organization of the Human Prostaglandin EP₃ Receptor Subtype Gene (PTGER3). *Genomics* 40: 425-434
3. Kunikata T *et al.* (2005) Suppression of allergic inflammation by the prostaglandin E receptor subtype EP₃. *Nat. Immunol.* 6: 524-531.
4. Narumiya S and FitzGerald GA (2001) Genetic and pharmacological analysis of prostanoid receptor function. *J. Clin. Invest.* 108: 25-30.
5. Ushikubi F *et al.* (1998) Impaired febrile response in mice lacking the prostaglandin E receptor subtype EP₃. *Nature* 395: 281-284.

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™ EP3 Prostanoid Receptor Frozen Cells, Product No. HTS092LRTA**

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.