

human Serotonin 5-HT_{1A} Receptor Cell Line

Product No.: ES-310-C

Lot No.: 1826798

Material Provided

Cells: 2 x 1 mL frozen aliquot (ES-310-CV)
Format: ~2.5 x 10⁶ cells /mL in freezing medium

Product Information

Cellular Background: CHO-K1

Cell Line Development: Our proprietary bicistronic expression plasmid containing the sequence coding for the human Serotonin 5-HT_{1A} receptor was transfected in CHO-K1 cells. Geneticin-resistant clones were obtained by limit dilution and compared for receptor expression levels using a radioligand binding assay. The clone with the highest receptor expression level was selected for characterization in binding and functional assays.

DNA Sequence: Identical to coding sequence of GenBank M83181.1. with the exception of a silent mutation in codon # 117 (GTG becomes GTC, both coding for a Val).

Corresponding Protein Sequence: Identical to GenBank NP_000515.2

Receptor expression level (B_{max}): Estimated to be 3.7 ± 2.3 pmol/mg protein, using [³H] 8-OH-DPAT

K_d for the above radioligand: 0.85 ± 0.49 nM

Shipping Conditions: Shipped on dry ice. Please ensure dry ice is still present in the package upon receipt or contact customer support.

Storage Conditions: Store in liquid nitrogen (vapor phase) immediately upon receipt.



Quality Control

The EC₅₀ for a reference agonist was determined in a LANCE *Ultra* cAMP assay performed on an EnVision® instrument. A mycoplasma test was performed using MycoAlert® (Lonza) mycoplasma detection kit. We certify that these results meet our quality release criteria.

5-Carboxamidotriptamine (5-CT) (EC₅₀): 0.74 nM

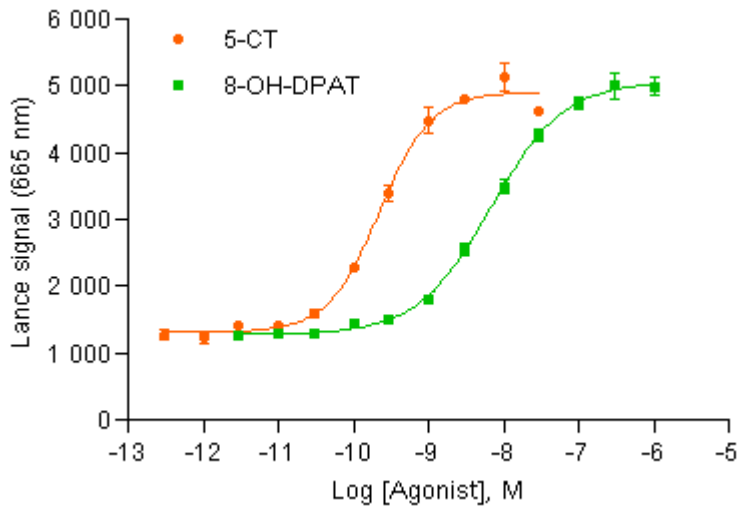
Stability: Cells were kept in continuous culture for at least 60 days and showed no decrease of receptor expression level in a saturation binding assay (stable B_{max} and K_d).

Mycoplasma: This cell line tested negative for mycoplasma.

Assay Procedures

We have shown for many of our GPCR cell lines that freshly thawed cells respond with the same pharmacology as cultured cells. All of our products validated in this way are available as frozen ready-to-use cells in our catalogue. This demonstrates that cells can be prepared and frozen in advance of a screening campaign simplifying assay logistics.

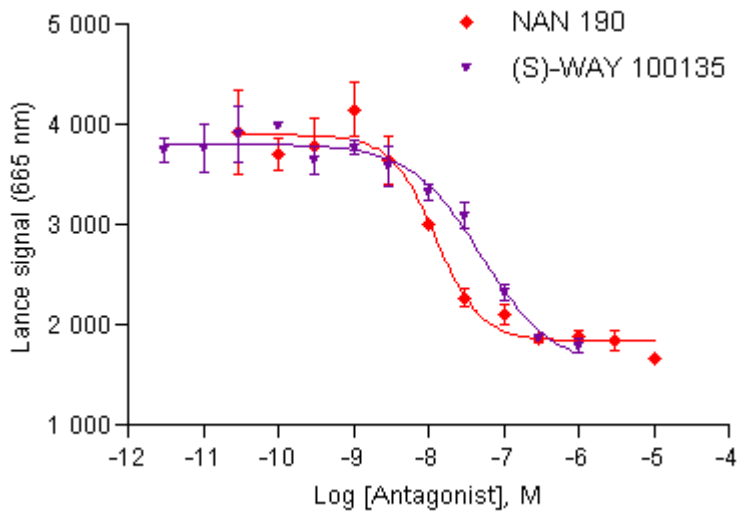
Typical Product Data -LANCE® Ultra cAMP Assay



Agonist	EC ₅₀ (M)
5-CT	2.2 x 10 ⁻¹⁰
8-OH-DPAT	6.7 x 10 ⁻⁹

Figure 1. Agonist Response in LANCE® Ultra cAMP assay

An agonist dose-response experiment was performed in 384-well format using 2500 cells/well. Frozen cells were thawed and incubated for 30-min with 10 µM Forskolin (FK) and the indicated agonist concentrations. Time-resolved fluorescence was measured on an EnVision® instrument. Data from a representative experiment are shown.



Antagonist	IC ₅₀ (M)
WAY100,135	1.3 x 10 ⁻⁸
NAN190	5.0 x 10 ⁻⁸

Figure 2. Antagonist Response in LANCE® Ultra cAMP assay

An antagonist dose-response experiment was performed in 384-well format using 2500 cells/well. Frozen cells were thawed and incubated for 30-min in the presence of 10 µM Forskolin (FK), a final concentration of 0.7 nM 5-CT, (corresponding to the EC₈₀), and the indicated antagonist concentrations. Time-resolved fluorescence was measured on an EnVision® instrument. Data from a representative experiment are shown.

Typical Product Data -Radioligand Binding Assay (Filtration)

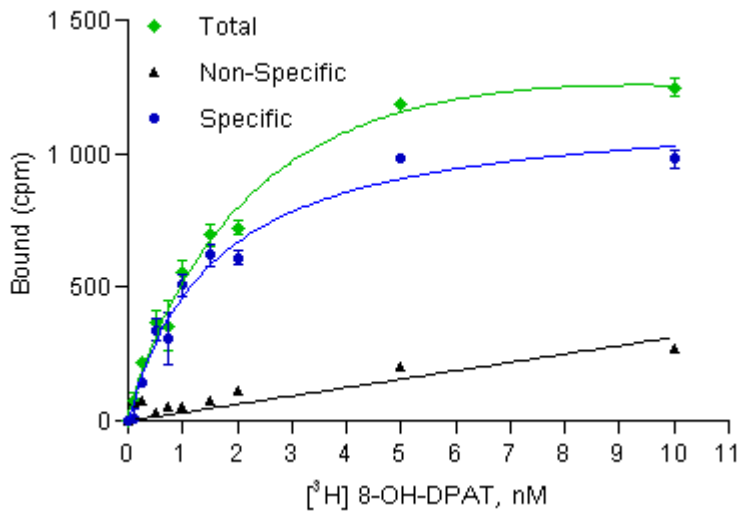
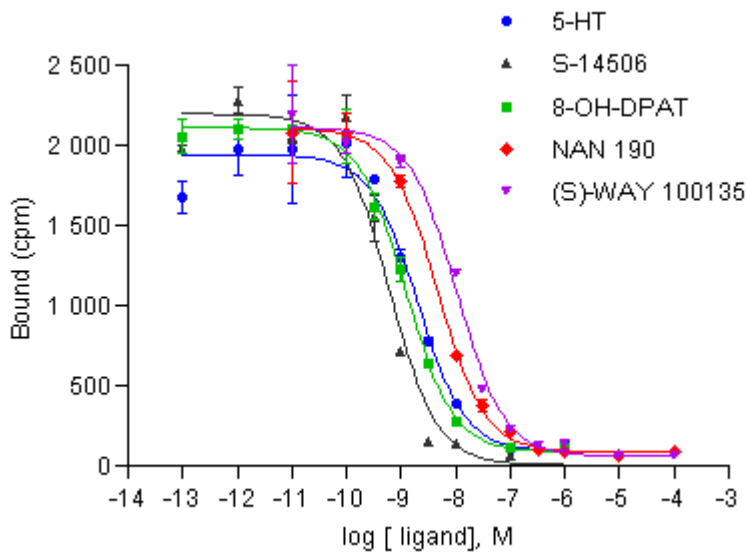


Figure 3: Saturation Binding Assay Curve (Filtration)

A saturation binding assay was performed in 96-well format using 2.5 µg membranes/well. Counts per minute (cpm) were measured on a TopCount® instrument. Data from a representative experiment are shown.

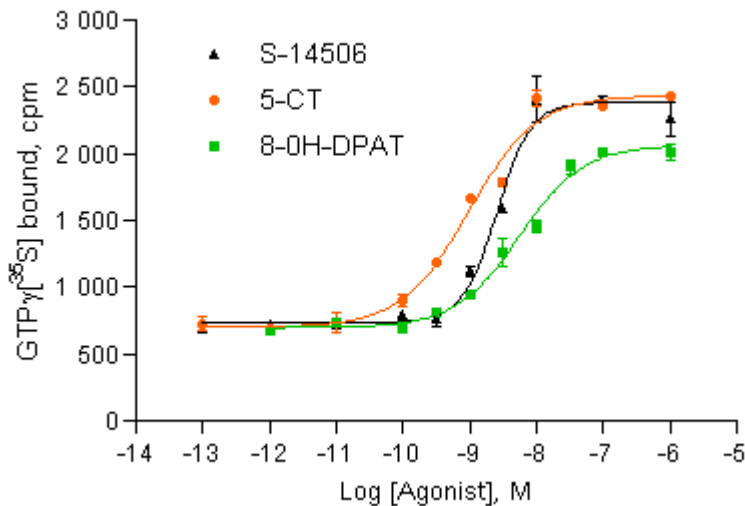


Agonist / Antagonist	IC ₅₀ (M)
5-HT	2.0 x 10⁻⁹
S-14506	6.1 x 10⁻¹⁰
8-OH-DPAT	1.2 x 10⁻⁹
NAN 190	4.6 x 10⁻⁹
(S)-WAY 100135	1.0 x 10⁻⁸

Figure 4: Competition Binding Assay Curve (Filtration)

A competition binding assay was performed in 96-well format using 2.5 µg membranes/well. Displacement of 1 nM [3H] 8-OH-DPAT was used. Counts per minute (cpm) were measured on a TopCount® instrument. Data from a representative experiment are shown.

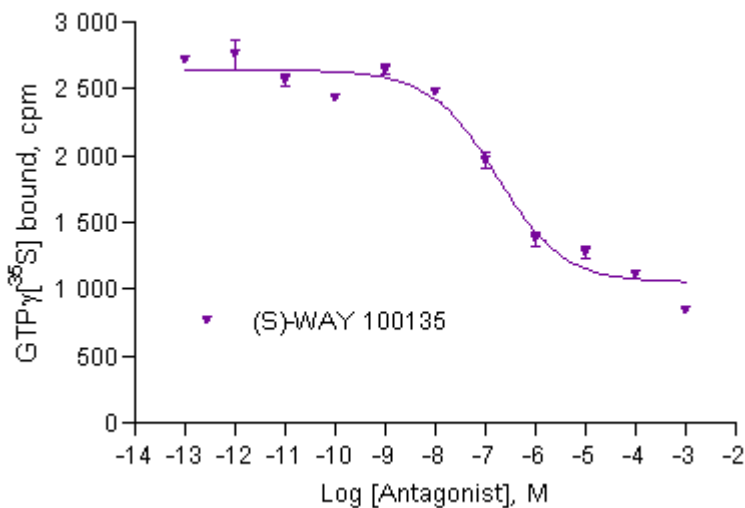
Typical Product Data - GTP γ S - SPA[®] Assay



Agonist	EC ₅₀ (M)
5-CT	9.7 x 10 ⁻¹⁰
8-OH-DPAT	5.5 x 10 ⁻⁹
S-14506	2.5 x 10 ⁻⁹

Figure 5. Agonist Response in GTP γ S - SPA[®] assay

An agonist dose-response scintillation proximity assay (SPA) was performed in 96-well format using 5 μ g membranes/well. Counts per minute (cpm) were measured on a TopCount[®] instrument. Data from a representative experiment are shown.



Antagonist	IC ₅₀ (M)
(S)-WAY 100135	1.7 x 10 ⁻⁷

Figure 6. Antagonist Response in GTP γ S - SPA[®] assay

An antagonist dose-response scintillation proximity assay (SPA) was performed in 96-well format using 5 μ g membranes/well. Counts per minute (cpm) were measured on a TopCount[®] instrument. Data from a representative experiment are shown.



LANCE® Ultra cAMP Assay Procedure

Stimulation Buffer: HBSS, 5 mM HEPES, 0.1 % Protease-free BSA, 0.5 mM IBMX, pH 7.4.

Cells/well: 2 500. For compounds not tested herein we recommend titrating the cells for optimal performance, i.e. 500-5 000 cells per assay point.

cAMP measurements were performed with the LANCE® Ultra cAMP 384 Kit (Revvity # TRF0262), according to the manufacturer instructions. Briefly:

Protocols for a 384-well white Optiplate (total assay volume of 20 µL):

cAMP Standard curve	G _s Agonist	G _s Antagonist	G _i Forskolin titration	G _i Agonist	G _i Antagonist
5 µL cAMP Standard	5 µL cell suspension	5 µL cell suspension	5 µL cell suspension	5 µL cell suspension	5 µL cell suspension
5 µL Stimulation Buffer	5 µL Agonist	2.5 µL Agonist	5 µL Forskolin	2.5 µL Forskolin	2.5 µL Forskolin/Agonist
-	-	2.5 µL Antagonist	-	2.5 µL Agonist	2.5 µL Antagonist
Incubate 30 min at room temperature (optional step for cAMP Standard curve)					
5 µL 4X Eu-cAMP Tracer Working Solution					
5 µL 4X ULight-anti-cAMP Working Solution					
Incubate 1 h at room temperature					
Read on an EnVision® instrument. Remove microplate seal prior to reading					

1. Cells in mid-log phase, grown in media without antibiotics for 18 hours prior to the experiment, were detached by gentle flushing with PBS-EDTA, recovered by centrifugation and resuspended in stimulation Assay Buffer at the concentration of 6.0×10^5 cells/mL.
2. Prepare the 4X Tracer Working Solution by making a 1/50 dilution of the Eu-cAMP stock solution in the cAMP Detection Buffer.
3. Prepare an ULight-anti-cAMP Intermediate Solution by making a 1/10 dilution of the ULight-anti-cAMP stock solution in cAMP Detection Buffer. Prepare the 4X ULight-anti-cAMP Working Solution by making a 1/30 dilution of the ULight-anti-cAMP intermediate solution in the cAMP Detection Buffer.

Notes:

For 96- and 1536-well formats, adjust proportionally the volume of each assay component in order to maintain the volume ratios for the 384-well format. Do not modify the Eu-cAMP and/or the ULight-anti-cAMP concentrations.



Membrane Radioligand Binding Assay Procedure (Filtration)

Note: The following are recommended assay conditions and may differ from the conditions used to generate the typical data shown in the above section.

Assay Buffer: 50 mM Tris-HCl pH 7.4, 10 mM MgSO₄, 0.5 mM EDTA, 0.1% ascorbic acid

Wash Buffer: Tris-HCl 50 mM pH 7.4 (ice cold)

Radioligand: [³H] LSD (Revvity # NET638)

Filters: Unifilter 96 GF/C (Revvity #6055690)

Membrane Binding Protocol:

Binding assays were performed in 550 µL total volume according to the following conditions. All dilutions are performed in assay buffer:

1. Membrane dilution:	2.5 µg of membranes per well, diluted in order to dispense 500 µL/well. Keep on ice.
2. Assembly on ice (in 96 Deep well plate)	<ul style="list-style-type: none">• 25 µL of assay buffer or of unlabeled ligand (Metergoline, 10 µM final) for determination of non specific binding• 25 µL of radioligand at increasing concentrations (see figure 3)• 500 µL of diluted membranes
Saturation Binding:	
Competition Binding:	<ul style="list-style-type: none">• 25 µL competitor ligand at increasing concentrations (see figure 4)• 25 µL of radioligand (1 nM final)• 500 µL of diluted membranes
3. Incubation:	60 min at 27°C.
4. Filters preparation:	GF/C filters were presoaked in 0.3% PEI at room temperature for at least 30 min.
5. Filtration:	Aspirate and wash 9 x 500 µL with ice cold wash buffer using a FilterMate Harvester.
6. Counting:	Add 30 µL/well of MicroScint™-O (Revvity # 6013611), cover filter with a TopSeal-A PLUS (Revvity # 6050185) and read on a TopCount®.



GTP γ S – SPA[®] Assay Procedure

Assay Buffer:	20 mM HEPES pH 7.4, 100 mM NaCl, 10 μ g/ml saponin, 3 mM MgCl ₂ ,
GDP concentration:	3 μ M GDP (final)
SPA Beads:	PVT-WGA (Revvity # RPNQ0001), 0.5 mg/well
Radioligand:	GTP γ S, [³⁵ S] - (Revvity # NEG030H)
Membranes:	5 μ g/well
Format:	96-well
Final volume:	100 μ L/well

GTP γ S-SPA assays were performed in 100 μ L total volume according to the following conditions. All dilutions are performed in assay buffer:

1. Membrane Dilution:	5 μ g of membranes per well, diluted in order to dispense 20 μ L/well. Keep on ice.
2. GDP preparation:	Prepare a 5-fold concentrated GDP solution (i.e. 15 μ M).
3. GTP γ S, [³⁵ S] - dilution:	Dilute GTP γ S, [³⁵ S] - to give ~25.000 dpm/20 μ L
4. Beads:	Dilute beads to 12.5 mg/mL (0.5 mg/20 μ L)
5. Assembly (in Optiplate™), Agonist Assay:	<ul style="list-style-type: none"> • 20 μL of 5x GDP dilution • 20 μL of 5x agonist dilutions at increasing concentrations • 20 μL of diluted membranes
Antagonist Assay:	<ul style="list-style-type: none"> • 20 μL of 5x GDP dilution • 20 μL of a 5x antagonist at increasing concentrations: 5x reference agonist dilution (to reach a final concentration corresponding to its EC₈₀) • 20 μL of diluted membranes
6. Pre-incubation:	Incubate for 15 min at room temperature (RT)
7. Assemble (continued)	<ul style="list-style-type: none"> • 20 μL of the GTPγS, [³⁵S] - dilution • 20 μL of the SPA Beads dilution
8. Incubation:	<ul style="list-style-type: none"> • Cover plate with a TopSeal, • Shake on an orbital shaker for 2 min, • Incubate for 30 min at RT • Centrifuge the plate for 10 min. at 2000 rpm, • Incubate for 0h to 1h at RT
9. Counting	Count for 1 min on a TopCount [®]

References

1. Kobilka B.K., Frielle T., Collins S., Yang-Feng T., Kobilka T.S., Francke U., Lefkowitz R.J. & Caron M.G. (1987) An intronless gene encoding a potential member of the family of receptors coupled to guanine nucleotide regulatory proteins. *Nature* 329:75-79.
2. Newman-Tancredi A., Wootton R. & Strange P.G. (1992) High-level stable expression of recombinant 5-HT_{1A} 5-hydroxytryptamine receptors in Chinese hamster ovary cells. *Biochem. J.* 285:933-938.
3. Raymond J.R., Albers F.J. & Middleton J.P. (1992) Functional expression of human 5-HT_{1A} receptors and differential coupling to second messengers in CHO cells. *Naunyn-Schiedeberg's Archs Pharmac.* 346:127-137.
4. Boess, F.G. and Martin, I.L. (1994). Molecular biology of 5-HT receptors. *Neuropharmacology* 33:275-317.
5. Xie D.-W., Deng Z.L., Ishigaki T., Nakamura Y., Suzuki Y., Miyasato K., Ohara K. & Ohara K. (1995) The gene encoding the 5-HT_{1A} receptor is intact in mood disorders. *Neuropsychopharmacology* 12:263-268.
6. Newman-Tancredi A. (2010) The importance of 5-HT_{1A} receptor agonism in antipsychotic drug action: rationale and perspectives. *Curr Opin Investig Drugs*

Materials and Instrumentation

The following tables provide the references of compounds and reagents used or recommended for the characterization of the human Serotonin 5-HT_{1A} receptor ValiScreen® cell line, as well as some advice on how to use these compounds:

Table 1. References of compounds used for functional characterization and binding assays

Name	Provider	Cat no	Working Stock Solution
5-Carboxamidotriptamine (5-CT)	Sigma	C117	10 mM in dH ₂ O
5-Hydroxytryptamine (5-HT)	Sigma	H-9523	10 mM in dH ₂ O
8-Hydroxy-DPAT (8-OH-DPAT)	Tocris	0529	10 mM in dH ₂ O
NAN 190	Sigma	N3529	10 mM in DMSO
(S)-WAY 100135	Tocris	1253	10 mM in DMSO
S-14506	Tocris	1771	10 mM in DMSO
[³ H] 8-OH-DPAT	Revvity	NET929	N/A
[³ H] LSD	Revvity	NET638	N/A

Table 2. References of cell culture media and assay buffers

Name	Provider	Cat no
HAM's F-12	Hyclone	SH30026.02
DMEM	Hyclone	SH30022.02
Advanced DMEM/F12 (serotonin receptors)	Invitrogen	12634-010
EMEM	BioWitthaker	06-174G
EX-CELL DHFR ⁻ media (DHFR deficient cell lines)	Sigma	C8862
FBS	Wisent	80150
FBS dialyzed	Wisent	80950
G418 (geneticin)	Wisent	400-130-IG
Zeocin	Invitrogen	R25005
Blasticidin	invitrogen	R210-01
Puromycin	Wisent	400-160-EM
Standard HBSS (with CaCl ₂ and MgCl ₂)	GIBCO	14025
HEPES	MP Biomedicals, LLC	101926
BSA, Protease-free	Sigma	A-3059
PEI	Sigma	P3143
Trypsin-EDTA	Hyclone	SH30236.02
Sodium Pyruvate	GIBCO	11360
L-Glutamine	GIBCO	25030
NEAA (non-essential amino acids)	GIBCO	11140
Forskolin	Sigma	F6886

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.