

Version: 02

Update (MM/DD/YYYY): 03/25/2025

Human Recombinant μ -Opioid Receptor Stable Cell Line

Cat. No. M00304

I. INTRODUCTION.....	1
II. BACKGROUND.....	2
III. REPRESENTATIVE DATA.....	2
IV. THAWING, SUBCULTURING AND CRYOPRESERVATION.....	4
V. REFERENCES.....	5
LIMITED USE LICENSE AGREEMENT.....	6

I. INTRODUCTION

Recombinant CHO-K1 cells stably overexpress human opioid receptor mu 1 (OPRM1) on the surface and contain high levels of G protein Gai and Gaq to couple with the receptor in downstream signaling pathways.

Catalog Number: M00304

Cell Line Name: CHO-K1/OPRM1/G α 15

Gene Synonyms: LMOR, M-OR-1, MOP, MOR, MOR1, OPRM

Expressed Gene: NCBI reference sequence NM_000914; no expressed tags

Target Protein: NP_000905

Host Cell: CHO-K1

Size: Two vials of frozen cells ($>1 \times 10^6$ per vial in 1 mL)

Culture Properties: Adherent

Freeze Medium: 45% Ham's F-12K (Kaighn's) (Cat. No. 21127, Life Technologies), 45% FBS (Cat. No. 10099-141, Life Technologies), 10% DMSO (Cat. No. D2650, Sigma)

Complete Growth Medium: Ham's F-12K (Kaighn's), 10% FBS

Culture Medium: Ham's F-12K (Kaighn's), 10% FBS, 200 μ g/ml Zeocin (Cat. No. R250-01, Life Technologies), 100 μ g/ml Hygromycin B (Cat. No. 10687010, Life Technologies)

Stability: Stable through more than 16 passages with no significant changes in assay performance or expression profile.

Application: Calcium flux assay, IP-One assay and cAMP accumulation assay.

Mycoplasma Status: Negative. The mycoplasma test was performed with MycoAlert™ PLUS Mycoplasma Detection Kit (Cat. No. LT07-318, Lonza).

Storage: Store cells in liquid nitrogen immediately upon receipt. Thaw and recover cells within one year from the date received.

II. BACKGROUND

The μ -opioid receptor (oprml) is the principal site of action in the brain by which morphine, other opiate drugs of abuse, and endogenous opioid peptides effect analgesia and alter mood. Opioid receptors belong to the rhodopsin family of G protein-coupled receptors (GPCRs). The three types of opioid receptors (μ , δ , and κ) have been shown to associate with each other in a homotypic or heterotypic fashion when expressed in heterologous cells. A member of the seven-transmembrane domain (TM) G protein-coupled receptor (GPCR) superfamily, the μ -opioid receptor modulates ion channels and second messenger effectors in an opioid agonist-dependent fashion that is reversible by the classic opiate antagonist naloxone.

III. REPRESENTATIVE DATA

Calcium Flux Analysis

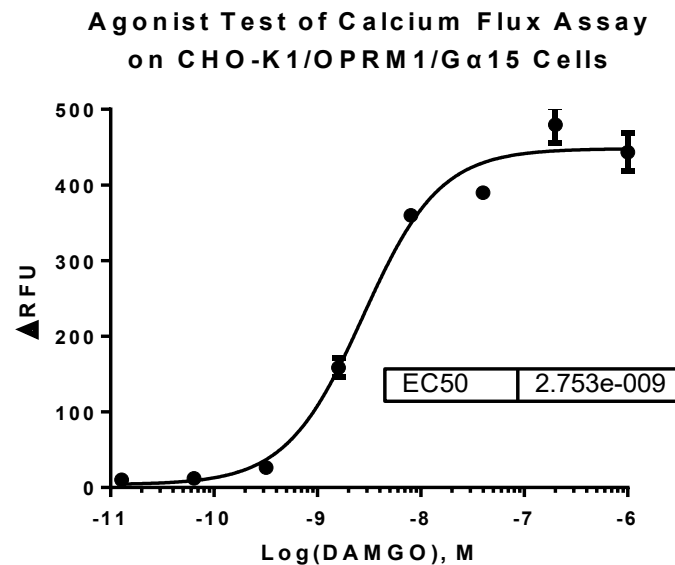


Figure 1. DAMGO-induced concentration-dependent stimulation of intracellular calcium mobilization in CHO-K1/OPRM1/G α 15 cells. The cells were loaded with Calcium-4 (Cat. No. R8142; Molecular Devices) prior to stimulation with OPRM1 agonist, DAMGO. The intracellular calcium change was measured by FLIPR^{TETRA}. The relative fluorescent units (RFU) were recorded and normalized to plot against the log of the cumulative doses of DAMGO (mean \pm SEM, n = 3). The EC₅₀ of DAMGO on CHO-K1/OPRM1/G α 15 cells was 2.75 nM.

Notes:

EC₅₀ value is calculated with four parameter logistic equation:

$$Y = \text{Bottom} + (\text{Top} - \text{Bottom}) / (1 + 10^{((\text{LogEC}_{50} - X) * \text{Hill Slope}))}$$

X is the logarithm of concentration. Y is the response.

Y is RFU and starts at Bottom and goes to Top along a sigmoid curve.

cAMP Accumulation Assay

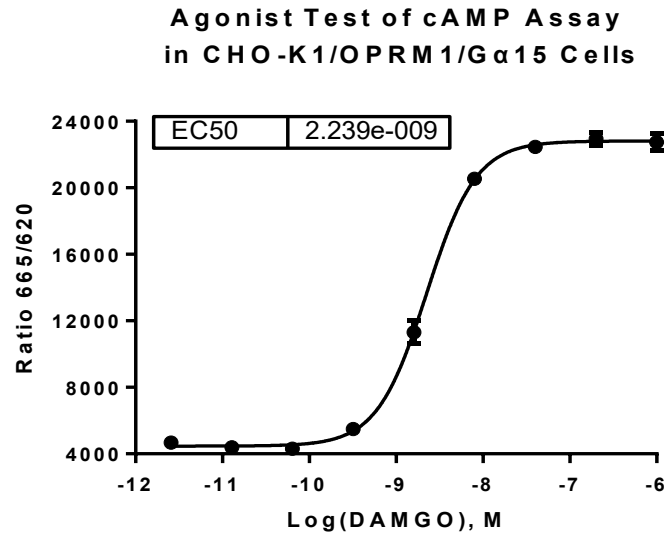


Figure 2. Dose dependent stimulation of intracellular cAMP accumulation upon treatment with DAMGO in CHO-K1/OPRM1/G α 15 cells. d2 acceptor fluorophore -labeled cAMP (Cat. No. 62AM4PEC; Revvity) and intracellular cAMP in CHO-K1/OPRM1/G α 15 cells competitively bind with Europium Cryptate-labeled anti-cAMP monoclonal antibody. The FRET signal decreases as the intracellular cAMP concentration rises and was measured by plate reader (Pherastar, BMG). The EC₅₀ of DAMGO on CHO-K1/OPRM1/G α 15 cells was 2.24 nM.

IP-One Accumulation Assay

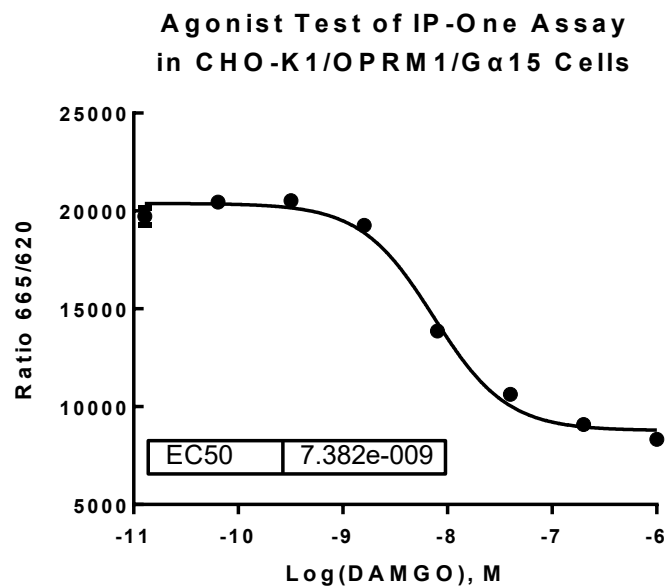


Figure 3. Dose dependent stimulation of intracellular IP-One accumulation upon treatment

with **DAMGO in CHO-K1/OPRM1/Gα15 cells**. d2 acceptor fluorophore-labeled IP-One (Cat. No. 62IPAPEB; Revvity) and intracellular IP-One in CHO-K1/OPRM1/Gα15 cells competitively bind with Europium Cryptate-labeled anti-IP-One antibody. The FRET signal decreases as the intracellular IP-One concentration rises and was measured by plate reader (Pherastar, BMG). The EC₅₀ of DAMGO on CHO-K1/OPRM1/Gα15 cells was 7.38 nM.

IV. THAWING AND SUBCULTURING

Thawing Protocol

1. Remove the vial containing the frozen cells from liquid nitrogen tank and place into a 37°C water bath immediately.
2. Thaw the cells quickly (within 1-2 minutes) by gently swirling the vial. Do not vortex the cells.
3. When the cells are almost completely thawed, take the vial out of the water bath and decontaminate it with 70% ethanol.
4. In a biosafety hood, transfer the cells to a sterile 15 ml conical tube. Add 9 ml of complete growth medium to the cells.
5. Pellet cells by centrifugation at 200 × g for 3-5 minutes at room temperature.
6. Carefully remove the supernatant with a pipette, leaving a small amount of medium to ensure that the cell pellet is not disturbed.
7. Resuspend the cells by gently flicking the tube. Gently add in 1 ml of complete growth medium.
8. Transfer the cell suspension into a 10 cm culture dish containing 10 ml of complete growth medium.
9. Grow the cells in an incubator at 37°C with 5% CO₂.
10. The cells will attach the dish in about 2-4 days. Check the status of the cells every day and don't disturb the cells till most cells attach well.
11. Change the medium with culture medium when cells grow well.

Sub-culturing Protocol

1. Remove the culture medium from the cells.
2. Wash cells with sterile PBS to remove all traces of serum which contains trypsin inhibitors.
3. Add 0.25% Trypsin/EDTA (Cat. No. 25200, Gibco) solution to the culture dish and observe the cells under an inverted microscope until the cell layer has dispersed (usually within 3-5 minutes).

Notes: To avoid cells clumping, do not agitate the cells by hitting or shaking the dish while waiting for the cells to detach. If cells are difficult to detach, place the dish in a 37°C incubator for about 2 minutes.

4. Add 6-8 ml of complete growth medium to the culture dish, aspirate the medium with cells by gentle pipetting and then add into a sterile falcon tube.
5. Centrifuge the cells at 200 x g for 5 minutes, and remove the medium.
6. Resuspend the cells in culture medium and add the cell suspension to a new culture dish.

7. Grow the cells in an incubator at 37°C with 5% CO₂.

Notes:

Subcultivation Ratio: 1:3 to 1:8.

Medium Renewal: Every 2 to 3 days.

Cryopreservation Protocol

1. Remove the cell culture medium, wash the cells with PBS once (optional), gently add enough trypsin to cover the cells and incubate for approximately 2 minutes in a 37°C incubator.
2. Resuspend in cell culture medium and transfer into a sterile 50 ml conical tube.
3. Count the viable cells using a hemocytometer. If preferred, also determine the cell viability. Cell viability should be at least 90% for good cryopreservation.
4. Centrifuge the cells at about 200 × g for 5 minutes at room temperature to pellet cells. Remove the supernatant gently without disturbing the cell pellet.
5. Resuspend cells by adding freezing medium to the tube to the required cell density (2-5 × 10⁶ cells/ml for best results).
6. Aliquot 1 ml each into cryogenic storage vials and secure the lids.
7. Transfer the vials into a cryo-freezing container at room temperature and put into a -80°C freezer. The temperature inside the cryo-freezing container should decrease steadily by 1°C/minute.
8. After approximately 24 hours, remove the vials from the cyro-freezing container and transfer into liquid nitrogen for long term storage.

V. REFERENCES

1. Ho MK, New DC, Wong YH. (2002) Co-expressions of different opioid receptor types differentially modulate their signaling via G (16). *Neurosignals*, 11 (2): 115-22.
2. Raynor K, Kong H, Chen Y, Yasuda K, Yu L, Bell GI, Reisine T. (1994) Pharmacological characterization of the cloned kappa-, delta-, and mu-opioid receptors. *Mol Pharmacol*, 45 (2): 330-4.
3. Yu VC, Eiger S, Duan DS, Lameh J, Sadée W. (1990) Regulation of cyclic AMP by the mu-opioid receptor in human neuroblastoma SH-SY5Y cells. *J Neurochem*, 55 (4): 1390-6.

Limited Use License Agreement

This is a legal agreement between you (Licensee) and GenScript USA Inc. governing use of GenScript's stable cell line products and protocols provided to licensee. By purchasing and using the stable cell line, the buyer agrees to comply with the following terms and conditions of this label license and recognizes and agrees to such restrictions:

- 1) The products are not transferable and will be used at the site where they were purchased. Transfer to another site owned by buyer will be permitted only upon written request by buyer followed by subsequent written approval by GenScript.
- 2) The purchaser cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party.
- 3) The products sold by GenScript are for laboratory and animal research purposes only. The products are not to be used on humans, for consumption, or for any unlawful uses.

GenScript USA Inc. will not assert against the buyer a claim of infringement of patents owned or controlled by GenScript USA Inc. and claiming this product based upon the manufacture, use or sale of a clinical diagnostic, therapeutic and vaccine, or prophylactic product developed in research by the buyer in which this product or its components has been employed, provided that neither this product nor any of its components was used in the manufacture of such product. For information on the use of this product for other purposes, contact Marketing Department, GenScript USA Inc., 860 Centennial Ave., Piscataway, NJ 08854, U.S.A. Phone: 1-732-885-9188.

For support contact us at:

400-025-8686-5810 or 5256 (CN)

1-732-885-9188 (USA)

65-3159-1898 (Asia Pacific)

31(0)-71-569-0120 (Europe)

81-3-6811-6572 (Japan)

02-6925-6649 (Korea)

44(1865)679988 (United Kingdom)

Web: www.genscript.com

For research use only. Not intended for human or animal clinical trials, therapeutic or diagnostic use.

Manufacturer: Nanjing GenScript Biotech Co., Ltd. No. 28 Yongxi Road, Jiangning District, Nanjing, Jiangsu, China