
Human Recombinant Cannabinoid Receptor 2 Stable Cell Line

Cat. No. M00433

Version 08072020

I. Introduction	1
II. Background	1
III. Assay Development	2
IV. Thawing and Subculturing	2
V. References	3
VI. Limited Use License Agreement.....	4

I. INTRODUCTION

Catalog Number: M00433

Cell Line Name: CHO-K1/CB2

Gene Synonyms: CNR2; CB2; CX5

Expressed Gene: Genbank Accession Number NM_001841; no expressed tags

Host Cell: CHO-K1

Culture Properties: Adherent

Quantity: Two vials of frozen cells (>1×10⁶ per vial)

Stability: More than 16 passages

Application: Functional assay for CB2 receptor (cAMP assay)

Freeze Medium: 45% culture medium, 45% FBS (Cat. #10099-141, Gibco), 10% DMSO (Cat. #D2650, Sigma)

Complete Growth Medium: Ham's F-12K (Kaighn's) (Cat. #21127, Life Technologies), 10% FBS

Culture Medium: Ham's F-12K (Kaighn's), 10% FBS, 400 µg/ml G418 (Cat. #10131-035, Gibco)

Mycoplasma Status: Negative*

Storage: Liquid nitrogen immediately upon receipt

II. BACKGROUND

The Cannabinoid receptor 2 (CNR2) is a G_{i/o}-coupled GPCR expressed in spleen, tonsils, bone marrow and peripheral blood leukocytes. There is one report that cannabinoid-induced inhibition of helper T cell activation is lost in macrophages obtained from CNR2 knockout mice. CNR2 is a potential therapeutic target in the treatment of various disease conditions, such as pain, multiple sclerosis, vascular disease, Parkinson's disease, and other central nerve system disorders.

* The mycoplasma test was performed with MycoAlert™ PLUS Mycoplasma Detection Kit of Lonza.

III. ASSAY DEVELOPMENT

This cell based assay is based on HTRF[®] technology (Homogeneous Time-Resolved Fluorescence). It is a competitive immunoassay that uses cAMP labeled with the d2 acceptor fluorophore and an anti-cAMP antibody labeled with Europium Cryptate. The FRET signal decreases as cAMP concentration rises.

Agonist Assay Protocol (According To Cisbio Document Reference: 62am4peb)

1. Seed 5 μ l CHO-K1/CB2 cells into a 384-well low volume plate, 3,000 cells per well.
2. Add 5 μ l compound (diluted in buffer with 0.25% DMSO) to each well and incubate the plate for 30 min at RT.
3. Add 5ul of cAMP-d2 conjugate solution to each well.
4. Add 5 μ l of cAMP-AB lysis buffer solution to each well.
5. Incubate the plate in the dark for one hour at RT.
6. Read the plate Varioskan (Thermo).

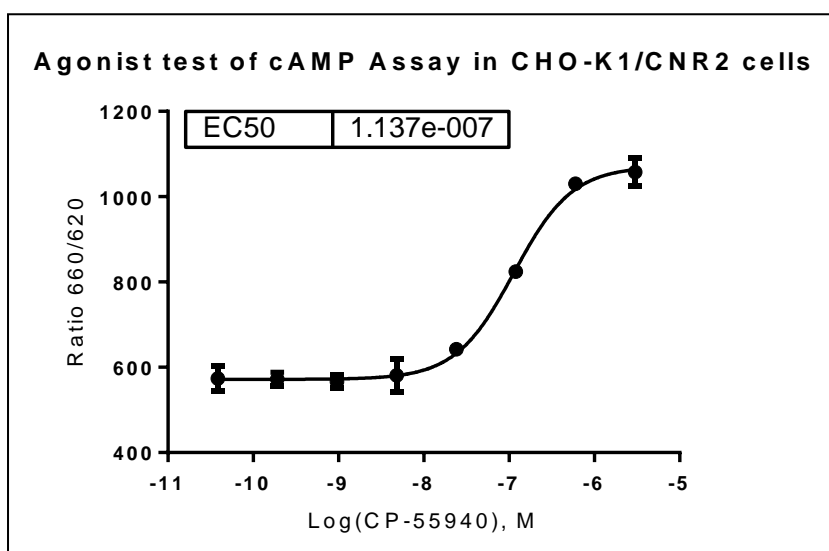


Figure 1. CP-55940-induced concentration-dependent stimulation of intracellular cAMP in CHO-K1/CNR2 cells. The EC₅₀ of CP-55940 on this cell was 0.11 μ M. Ratio = A660nm/B620nm x 10⁴

IV. THAWING AND SUBCULTURING

Thawing Protocol

1. Remove the vial from liquid nitrogen tank and thaw cells quickly in a 37°C water-bath.
2. Just before the cells are completely thawed, decontaminate the outside of the vial with 70% ethanol and transfer the cells to a 15 ml centrifuge tube containing 9 ml of complete growth medium.
3. Pellet cells by centrifugation at 200 x g force for 5 min, and discard the medium.
4. Resuspend the cells in complete growth medium.
5. Add 10 ml of the cell suspension in a 10 cm dish.
6. Add G418 to concentrations of 400 μ g/ml the following day.

Sub-culturing Protocol

1. Remove and discard culture medium.
2. Wash cells with PBS (pH=7.4) to remove all traces of serum that contains trypsin inhibitor.
3. Add 2.0 ml of 0.05% (w/v) Trypsin- EDTA (GIBCO, Cat No. 25300) solution to 10 cm dish and observe the cells under an inverted microscope until cell layer is dispersed (usually within 3 to 5 minutes).

Note: To avoid clumping, do not agitate the cells by hitting or shaking the dish while waiting for the cells to detach. Cells that are difficult to detach may be placed at 37°C to facilitate dispersal.

4. Add 6.0 to 8.0 ml of complete growth medium and aspirate cells by gently pipetting, centrifuge the cells 200 x g force for 5min, and discard the medium.
5. Resuspend the cells in culture medium and add appropriate aliquots of the cell suspension to new culture vessels.
6. Incubate cultures at 37°C.

Subcultivation Ratio: 1:3 to 1:8

Medium Renewal: Every 2 to 3 days

V. REFERENCES

1. T Howlett AC et al. International Union of Pharmacology. XXVII. Classification of cannabinoid receptors. Pharmacol Rev. 2002 Jun; 54(2):161-202.
2. Buckley NE et al. Immunomodulation by cannabinoids is absent in mice deficient for the cannabinoid CB (2) receptor. Eur J Pharmacol. 2000 May 19; 396(2-3):141-9.

GenScript USA Inc,
860 Centennial Ave.
Piscataway, NJ 08854
Toll-Free: 1-877-436-7274
Tel: 1-732-885-9188, Fax: 1-732-210-0262
Email: product@genscript.com
Web: <http://www.genscript.com>
For Research Use Only.

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. Fax: 1-732-210-0262. Email: marketing@genscript.com.