

## MULTISCREEN™ STABLE CELL LINE MOUSE RECOMBINANT CB1 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** Cm1229-1A

**Lot Number:** Cm1229-1A-112316

**Quantity:** 1 vial ( $2 \times 10^6$ ) frozen cells

**Freeze Medium:** 90% FBS, 10% DMSO

**Host cell:** CHO-K1

**Transfection:** Expression vector containing full-length *Mus musculus* CB1 cDNA (GenBank Accession Number: BC079564) with FLAG tag sequence at N-terminus.

**Recommended Storage:** Liquid nitrogen upon receiving

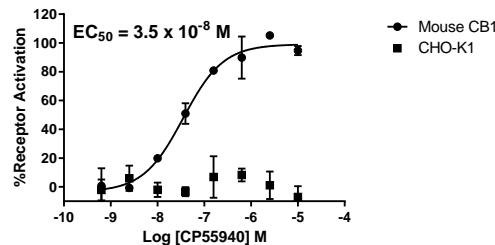
**Propagation Medium:** DMEM/F12, 10% FBS, 10  $\mu$ g/mL puromycin

**Stability:** In progress

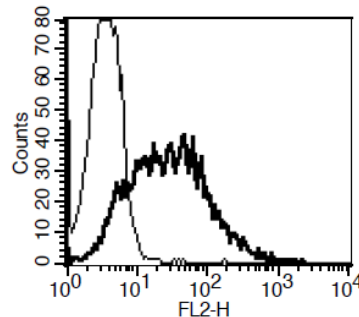
**Background:** Cannabinoid Receptor 1, CNR1 also known as CB1, is involved in cannabinoid induced CNS effects. It acts by inhibiting intracellular adenylate cyclase activity and could be a receptor for anandamide. CNR1 is a potential target for the development of novel therapeutic drugs in the treatment of various conditions, such as pain, feeding disorders, vascular disease, Parkinson's disease, and other central nerve system disorders.

**Application:** Functional assays

**Figure 1:**



**Figure 2:**



**Figure 1:** Dose-dependent stimulation of intracellular cAMP level upon treatment with ligand., measured with Multiscreen™ TR-FRET cAMP Assay Kit (Multispan MSCM01). **Figure 2:** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

#### References:

Gerard, C., C. Mollereau, et al. (1990). "Nucleotide sequence of a human cannabinoid receptor cDNA." *Nucleic Acids Res* 18(23): 7142

Mendizabal, V. E. and E. Adler-Graschinsky (2003). "Cannabinoid system as a potential target for drug development in the treatment of cardiovascular disease." *Curr Vasc Pharmacol* 1 (3): 301-13

**FOR RESEARCH USE ONLY.**

All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.