

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT CB1 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1229

**Lot Number:** C1229-062308

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

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human CNR1 cDNA (GenBank Accession Number NM\_016083.3) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

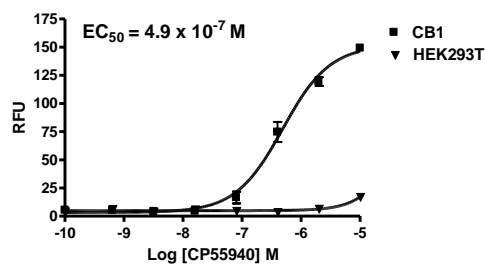
**Propagation Medium:** DMEM, 10% FBS, 1  $\mu\text{g}/\text{mL}$  puromycin

**Stability:** In progress

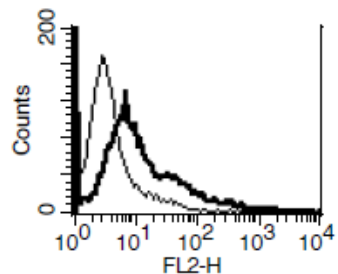
**Background:** Cannabinoid Receptor 1, CNR1 also known as CB1, is involved in cannabinoid induced CNS effects. It acts by inhibiting intracellular adenylate cyclase activity and could be a receptor for anandamide. CNR1 is a potential target for the development of novel therapeutic drugs in the treatment 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 calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **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:

Mendizabal and Adler-Graschinsky (2003) Cannabinoid system as a potential target for drug development in the treatment of cardiovascular disease. *Curr Vasc Pharmacol* 1:301-313.

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

**FOR RESEARCH USE ONLY.**

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