

**MULTISCREEN™ DIVISION ARRESTED CELL LINE  
HUMAN RECOMBINANT CB1 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1229-1a

**Lot Number:** DC1229-1a-021012

**Quantity:** 1 vial ( $4 \times 10^6$ ) frozen cells

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

**Host cell:** CHO-K1

**Transfection:** Expression vector containing full-length human CB1 cDNA (GenBank Accession Number: NM\_016083) with FLAG tag sequence at N-terminus.

**Recommended Storage:** Liquid nitrogen upon receiving

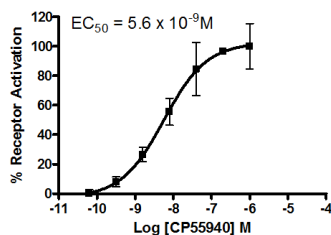
**Propagation Medium:** DMEM/F12, 10% FBS

**Stability:** Stable for 1-2 days after thawing

**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 1.** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP accumulation upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

**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.**

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