

MULTISCREEN™ STABLE CELL LINE MOUSE RECOMBINANT CB1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: Cm1229-1B

Lot Number: Cm1229-1B-112316

Quantity: 1 vial (2×10^6) 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 μ 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

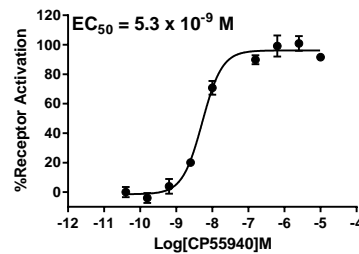


Figure legend: Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP Assay Kit (Multispan MSCM01)

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.

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