

MULTISCREEN™ DIVISION-ARRESTED CELL LINE HUMAN RECOMBINANT CB2 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1230-1a

Lot Number: DC1230-1a-021925

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Cellbanker 2

Host cell: CHO-K1

Transfection: Expression vector containing full-length human CNR2 cDNA (GenBank Accession Number NM_001841) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS

Background: Cannabinoid receptor 2 (CB2 or CNR2) is involved in cannabinoid-induced central nerve system effects. It could be a receptor for anandamide. The receptor 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.

Application: Functional assays

Figure 1

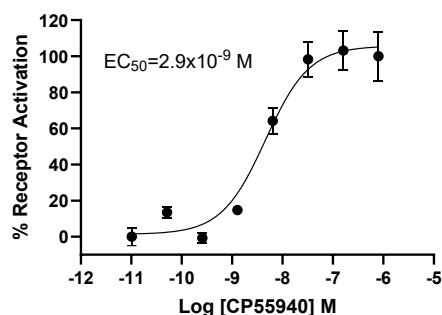


Figure 1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with MULTISCREEN™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

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.