

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT CB1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: MC1229

Lot Number: MC1229-100721

Quantity: 1 vial (8.15mg/mL),
1, 5, 10mg

Packaging Buffer: 20mM Gly-Gly, 1
mM MgCl₂, 25mM Sucrose (pH 7.2)

Host cell: HEK293T

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 Avoid repeated
freeze-thaw.

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.

Figure 1

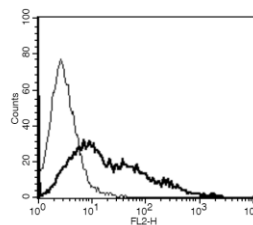


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

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.