

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT CB2 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1230-1a

Lot Number: C1230-1a-041421

Quantity: 1 vial (7.6mg/mL),
1 mg

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

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. Avoid
repeated freeze-thaw

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.

Figure 1

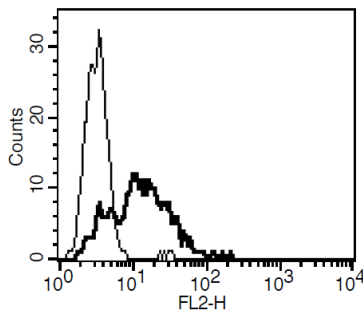


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

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.

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