

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT D4 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: MC1338

Lot Number: MC1338-020726

Quantity: Available 1mg, 5mg, or 10mg per vial

Protein concentration: 12.75 mg/mL

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

Host cell: HEK293T

Transfection: Full-length Human DRD4 cDNA (GenBank Accession NM_000797) with FLAG-tag sequence at the N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Background: The human dopamine receptor D4 is a D2-like receptor that inhibits adenylyl cyclase activity and activates K⁺ channels. D4 receptor antagonists show great potential in the treatment of human personality and psychiatric disorders, such as ADHD (Attention deficit hyperactivity disorder), schizophrenia, alcoholism, and drug addiction.

Application: Functional assays

Figure 1

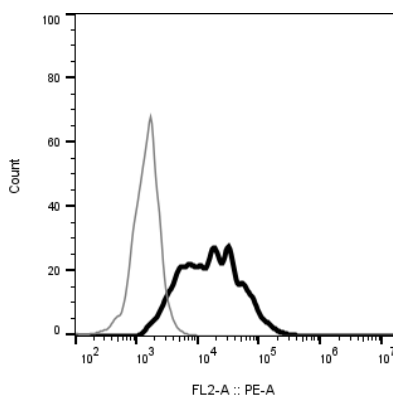


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Kulkarni SK, et al. (2000) Dopamine D4 receptors and development of newer antipsychotic drugs. *Fundam Clin Pharmacol* 14(6):529-39.

Missale C, et al. (1998) Dopamine receptors: from structure to function. *Physiol Rev* 78(1):189-225.

Van Tol HH, et al. (1991) Cloning of the gene for a human dopamine D4 receptor with high affinity for the antipsychotic clozapine. *Nature* 350(6319):610-4

FOR RESEARCH USE ONLY.

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