

**MULTISCREEN™ MEMBRANE PREPARATION
MOUSE RECOMBINANT mGLUR4 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: MHGm1191-1a

Lot Number: MHGm1191-1a-052215

Quantity: 7.07 mg/ml, 5mg

Packaging Buffer: 20 mM Gly-Gly, 1 mM MgCl₂, 250 mM Sucrose

Host cell: CHO Gαq5

Transfection: Expression vector containing full-length mouse GRM4 cDNA (GenBankAccession Number BC072635) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Background: L-glutamate is the major excitatory neurotransmitter in the central nervous system and activates both ionotropic and metabotropic glutamate receptors. The metabotropic glutamate receptors, which are G protein-coupled receptors, have been divided into 3 groups on the basis of sequence homology, putative signal transduction mechanisms and pharmacologic properties. Group II and group III receptors are linked to the inhibition of the cyclic AMP cascade, but differ in their agonist selectivity. Group III agonists include L-2-amino-4-phosphonobutyrate (L-AP4) and L-serine-O-phosphate.

Application: Functional assays

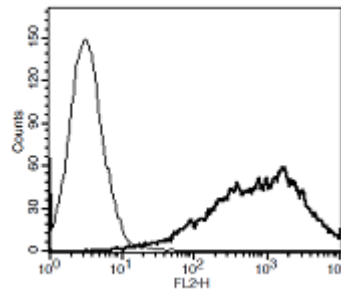


Figure 1. Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Wu *et al.* (1998) Group III human metabotropic glutamate receptors 4, 7 and 8: molecular cloning, functional expression, and comparison of pharmacological properties in RGT cells. *Mol Brain Res* 53:88-97.

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.