

PRODUCT DATASHEET

ChemiScreen™ Recombinant Human mGlu₂ Metabotropic Glutamate Receptor Membrane Preparation

CATALOG NUMBER:	HTS146M	QUANTITY:	200 units
LOT NUMBER:	21K1202	VOLUME/CONCENTRATION	1 mL, 1 mg/mL

BACKGROUND: Glutamate is a main excitatory neurotransmitter in the central nervous system, and it plays a role in learning, memory and neurotoxicity. The biological actions of glutamate are mediated by ionotropic and metabotropic glutamate receptors, which are ion channels and GPCRs respectively. Metabotropic glutamate receptors (mGluRs) are members of the class 3 G-protein coupled receptor family, which are characterized by a large extracellular domain. They are further classified into group I, II, and III mGluRs on the basis of their sequence identity, pharmacology, and signal transduction mechanism. Group I (mGlu₁ and mGlu₅) couple to the phospholipase C pathway through G_{αq}, whereas group II (mGlu₂ and mGlu₃) and group III (mGlu₄, mGlu₆, mGlu₇, and mGlu₈) negatively couple to the adenylyl cyclase pathway through G_{αi} (Conn and Pin, 1997). Agonists of the Group II metabotropic glutamate receptors, mGlu₂ and mGlu₃, display efficacy in animal models of anxiety and psychosis. A key role for mGlu₂ in mediating these effects is indicated by the observation that selective allosteric potentiator of mGlu₂ also retains antipsychotic-like activities in mice (Galici *et al.*, 2005). In addition, mGlu_{2/3} agonists display analgesic activity in animal models (Jones *et al.*, 2005). mGlu₂ membrane preparations are crude membrane preparations made from our proprietary stable recombinant cell lines to ensure high-level of GPCR surface expression; thus, they are ideal HTS tools for screening of mGlu₂ interactions with its ligands.

APPLICATIONS: Radioligand binding assay

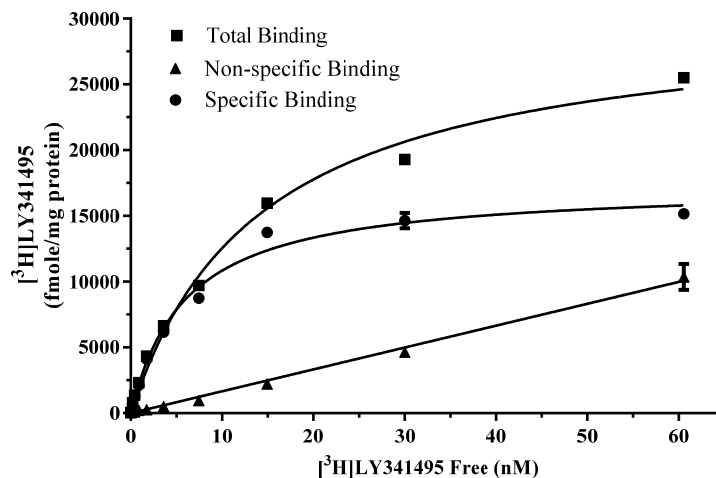


Figure 1. Saturation binding for mGlu₂ membrane preparation. 2 µg/well mGlu₂ Membrane Preparation was incubated with increasing amount of 3H-labeled LY341495 in the absence (total binding, TB) or presence (nonspecific binding, NSB) of unlabeled LY-354740 at 10µM. Specific binding (SB) was determined by subtracting NSB from TB.

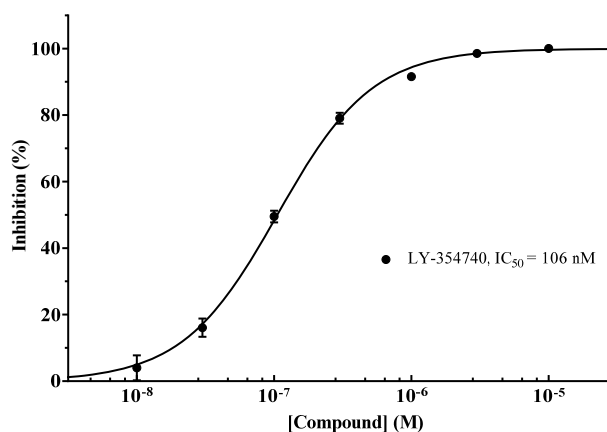


Figure 2. Competition binding for mGlu₂ membrane preparation. 2 µg/well mGlu₂ Membrane Preparation and wild-type Chem-1 Membrane Preparation (Chemicon catalog # HTS000MC1) were incubated in a 96-well plate with 2 nM ³H-labeled LY341495 and increasing concentrations of unlabeled LY-354740. More than 10-fold signal:background was obtained with LY-354740.

SPECIFICATIONS: 1 unit = 5 µg
 B_{max} for [³H] LY341495 binding: 17.4 pmol/mg protein
 K_d for [³H] LY341495 binding: ~6.0 nM
 Signal:background: >10-fold

TRANSFECTION: Human GRM2 cDNA encoding mGlu₂ (Accession number NM_000839)

HOST CELLS: Chem-1, an adherent cell line expressing the promiscuous G-protein, Gα15 and with no endogenous expression of mGlu₂.

RECOMMENDED ASSAY CONDITIONS: Membranes are mixed with radioactive ligand and unlabeled competitor (see Figures 1 and 2 for concentrations tested) in binding buffer in a nonbinding 96-well plate, and incubated for 1h. Prior to filtration, an GF/B 96-well harvest plate is coated with 10 mM KH₂PO₄, 100 mM KBr, pH 7.6. Binding reaction is transferred to the filter plate, and washed 4 times (~1 mL per well per wash) with Wash Buffer. The plate is dried and counted.

Binding buffer: 10 mM KH₂PO₄, 100 mM KBr, pH 7.6

Radioligand: [³H]-LY341495 (ARC, ART1439)

Wash Buffer: 10 mM KH₂PO₄, 100 mM KBr, pH 7.6

One package contains enough membranes for at least 200 assays (units), where a unit is the amount of membrane that will yield greater than 10-fold signal:background with ³H-labeled LY341495.

PRESENTATION: Liquid in packaging buffer: 50 mM Tris pH 7.4, 10% glycerol and 1% BSA with no preservatives.
 Packaging method: Membrane proteins were adjusted to the indicated concentration in 1 ml packaging buffer, rapidly frozen, and stored at -80°C.

STORAGE/HANDLING: Store at -70°C. Product is stable for at least 6 months from the date of receipt when stored as directed. Do not freeze and thaw.

REFERENCES:

1. Conn PJ and Pin JP (1997) Pharmacology and functions of metabotropic glutamate receptors. *Annu. Rev. Pharmacol. Toxicol.* 37: 205-37
2. Galici R *et al.* (2005) A selective allosteric potentiator of metabotropic glutamate (mGlu) 2 receptors has effects similar to an orthosteric mGlu2/3 receptor agonist in mouse models predictive of antipsychotic activity. *J. Pharmacol. Exp. Ther.* 315(3):1181-7
3. Jones CK *et al.* (2005) Analgesic effects of the selective group II (mGlu2/3) metabotropic glutamate receptor agonists LY379268 and LY389795 in persistent and inflammatory pain models after acute and repeated dosing. *Neuropharmacology* 49 Suppl 1:206-18.

FOR RESEARCH USE ONLY; NOT FOR USE IN DIAGNOSTIC PROCEDURES. NOT FOR HUMAN OR ANIMAL CONSUMPTION

Unless otherwise stated in our catalog or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals. No part of these works may be reproduced in any form without permission in writing.

Eurofins Pharma Bioanalytics Services US Inc. is an independent member of Eurofins Discovery Services