

**MULTISCREEN™ DIVISION ARRESTED CELL LINE  
HUMAN RECOMBINANT MGLUR5 RECEPTOR**

**PRODUCT INFORMATION**

**Catalog Number:** DC1192

**Lot Number:** DC1192-121610

**Quantity:** 1 vial ( $4 \times 10^6$ ) frozen cells

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human GRM5 cDNA (NM\_000842.1) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM with GlutaMAX (Gibco 10566), 10% FBS (dialyzed), 2 mM sodium pyruvate

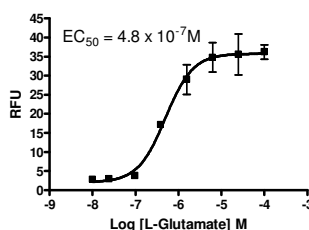
**Stability:** Stable for 2-3 days after thawing

**Data sheet**

**Background:** The metabotropic glutamate receptors (mGluRs) are divided into 3 groups on the basis of sequence homology, putative signal transduction mechanisms, and pharmacological properties. Group II and group III mGluRs are linked to inhibition of the cyclic AMP cascade, but differ in their agonist selectivity. mGluR5 is a member of Group I. It mainly expresses the brain although its precise role in various cortical functions remains unclear.

**Application:** Functional assays

**Figure 1**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation.

**References:**

Romano *et al.* (1995) Distribution of metabotropic glutamate receptor mGluR5 immunoreactivity in rat brain. *J Comp Neurol* 355:455-469.

She *et al.* (2009) Roles of mGluR5 in synaptic function and plasticity of the mouse thalamocortical pathway. *Eur J Neurosci* 29:1379-1396.

**FOR RESEARCH USE ONLY.**

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com  
sales@multispaninc.com  
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817  
Fax: +1 (510) 887-0863  
26219 Eden Landing Road  
Hayward, CA 94545-3718  
U.S.A.