

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT GAL1 RECEPTOR

PRODUCT INFORMATION

Catalog Number: CG1178-1

Lot Number: CG1178-1-021606

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1 Gαq5

Transfection: Expression vector containing full-length human GAL1 cDNA (GenBank Accession Number NM_001480) with flag tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM-F12, 10% FBS, 10 μg/mL puromycin, 150 μg/mL hygromycin

Stability: Stable after minimum of two months continuous growth

Data sheet

Background: The diverse physiological effects of Galanin, a biologically active neuropeptide, are mediated through cell surface G protein-coupled receptors. To date, three galanin receptor subtypes, GAL1, GAL2 and GAL3, have been cloned. Galanin, widely distributed in the central and peripheral nervous systems and the endocrine systems, binds to galanin receptors to induce several regulatory functions in neuronal cells, including neuroregeneration, control of endocrine and exocrine secretions, and modulation of sensory and behavioral functions. Galanin agonists have been shown to have therapeutic application in treatment of chronic pain; galanin antagonists have therapeutic potential in treatment of Alzheimer's disease, depression, and feeding disorders.

Application: Functional assays

Figure 1

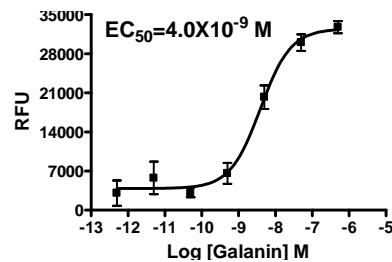


Figure 2

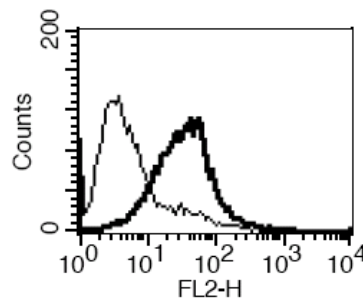


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Branchek *et al.* (1998) Molecular biology and pharmacology of galanin receptors. *Ann N Y Acad Sci* 863: 94-107.

Wang *et al.* (1998) Differential intracellular signaling of the GalR1 and GalR2 galanin receptor subtypes. *Biochemistry* 37:6711-6717.

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.