

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT GLP-2 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1268

Lot Number: C1268-122011

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human GLP-2 cDNA (GenBank Accession Number NM_004246) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μ g/mL puromycin

Stability: Stable in culture for minimum of two months

Background: Glucagon-like peptide 2 receptor belongs to the glucagon-secretin receptor superfamily of GPCRs. The human GLP-2 receptor gene is localized on chromosome 17p13.3. The GLP-2 receptor exhibits ~50% amino acid identity with GLP-1 receptor. GLP-2 receptor binds to Glucagon-like peptide-2 (GLP-2) is a nutrient-responsive hormone and activates adenylate cyclase pathway and to a lesser extent activates MAP kinases. GLP-2 receptors are found in the central nervous system and gastrointestinal tract, with the highest expression levels in jejunum. The principal role of GLP-2 receptors appears to be the maintenance of the growth, nutrient absorption, cell proliferation, apoptosis, mucosal blood flow and suppressing gastric motility and secretion. The regenerative and cytoprotective properties of GLP-2 contribute to its therapeutic potential for the treatment of patients with intestinal disease. Recent studies have suggested that GLP-2 not only modulates intestinal stem cell behavior but may also promote carcinogenesis in models of sporadic colon cancer. Further consideration of the molecular cross-talk and downstream signaling pathways mediating the intestinotropic effects of GLP-2 is important. A detailed delineation of the signaling pathways activated by gut peptide GPCRs, as exemplified by GLP-1 and GLP-2, may provide new therapeutic targets for the treatment of human disorders such as diabetes and intestinal disease, respectively

Application: Functional assays

Figure 1

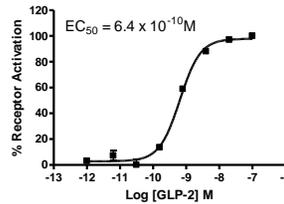


Figure 1. Dose-dependent stimulation of intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

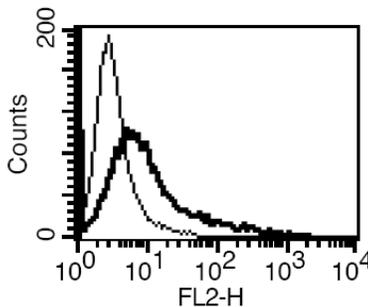
References:

Jennifer L. Estall and Daniel J. Drucker (2003) Dual Regulation of Cell Proliferation and Survival via Activation of Glucagon-Like Peptide-2 Receptor Signaling. *J. Nutr.* 133: 3708–3711.

Rocha *et al* (2004) Glucagon-like peptide-2: divergent signalling pathways. *J.Surg.Res.* 121: 5-9.

Brubaker and Drucker (2004) Glucagon-like peptides regulate cell proliferation and apoptosis in the pancreas, gut, and central nervous system. *Endocrinol.* 145: 2653-2659.

Figure 2



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.