

MULTISCREEN™ DIVISION ARRESTED CELL LINE HUMAN RECOMBINANT GPR37L1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1098

Lot Number: DC1098-042321

Quantity: 1 vial (4 x 10⁶) frozen cells

Freeze Medium: Cellbanker 2
(Amsbio)

Host cell: HEK293T

Transfection: Expression vector containing full-length human GPR37L1 cDNA (GenBank Accession Number NM_004767) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Background: GPR37L1 (G-protein coupled receptor 37 like 1) is also known as ETBR-LP-2 (Endothelin B receptor-like protein-2) and is an orphan receptor. The receptor is strongly expressed in the human central nervous system, e.g. in cerebellar Bergmann glia, cerebral cortex, and internal capsule fibers. The fragment of this receptor is detectable in human cerebrospinal fluid.

Application: Functional Analysis

Figure 1

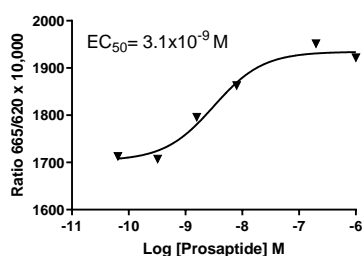


Figure 1. Dose-dependent inhibition of forskolin stimulated intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

References:

Valdenaire *et al.* (1998) A new family of orphan G protein-coupled receptors predominantly expressed in the brain. *FEBS Lett* 424:193-196.

Stark *et al.* (2001) Peptide repertoire of human cerebrospinal fluid: novel proteolytic fragments of neuroendocrine proteins. *J Chromatogr B Biomed Sci Appl* 754:357-367.

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.