

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT RXFP3 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1306

Lot Number: C1306-091609

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 RXFP3 cDNA (GenBank accession number: NM_016568.2) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: In progress

Background: The relaxin family peptide receptor RXFP3, also known as GPCR135, RLN3R1 or SALPR, is a 469-amino acid, 7-transmembrane protein. RT-PCR detected RXFP3 expression in various sections of the brain and the testis. RXFP3 is known to be important to several functions of the central nervous system, especially sensory processing under stressful conditions.

Application: Functional assays
Figure 1

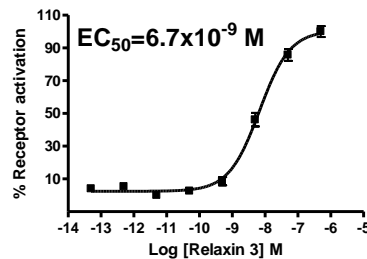


Figure 2

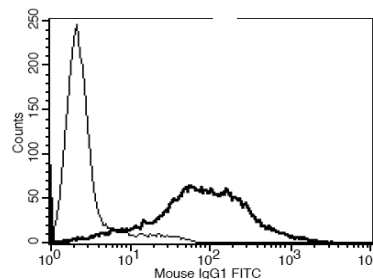


Figure 1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody.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

References:

Cgen *et al.* (2005) Pharmacological characterization of relaxin-3/INSL7 receptors GPCR135 and GPCR142 from different mammalian species. *J Pharmacol Exp Ther* 312:83-95.

Liu *et al.* (2003) Identification of relaxin-3/INSL7 as an endogenous ligand for orphan G-protein-coupled receptor GPCR135. *J Biol Chem* 278:50754-50764.

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.