

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT RXFP2 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1289NTb

Lot Number: C1289NTb-091015

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 RXFP2 cDNA (GenBank accession number: NM_021634)

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: Stable for minimum of two months in continuous growth

Background: The relaxin family peptide receptor RXFP2, also known as LGR8, is a 754-amino acid 7 transmembrane protein. RT-PCR detected RXFP2 expression in various organs of the reproductive system along with the brain, kidney, thyroid, muscle, peripheral blood cell, and bone marrow. RXFP2 is known to be important in supporting germ cell function in the testis and ovaries. Additionally, RXFP2 mutations are linked to osteoporosis in men, suggesting that it plays a central role in bone physiology.

Application: Functional assays

Figure 1

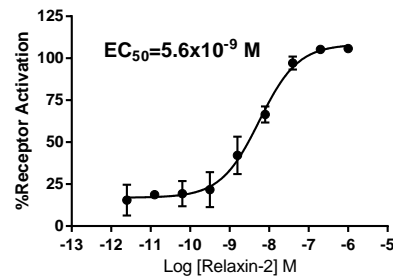


Figure 2

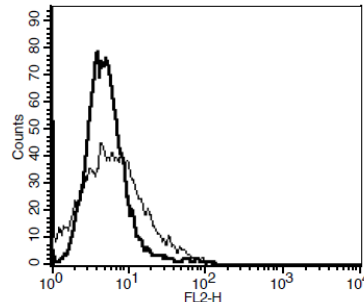


Figure 1. Dose-dependent increase of intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-RXFP2 antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Hall *et al.* (2006) Relaxin family peptide receptors, RXFP1 and RXFP2, modulate cAMP signalling by distinct mechanisms.. *Mol Pharmacol* (1):214-26.

Kumagi *et al.* (2002) INSL3/Leydig insulin-like peptide activates the LGR8 receptor important in testis descent.. *J Biol Chem* 277: 31283-31286.

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.