

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT RXFP2 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: MC1289NTa

Lot Number: MC1289NTa-020316

Quantity: 1 vial (10.5mg/mL, 1mg)

Packaging Buffer: 20 mM Tris-HCl (pH7.5), 2 mM EDTA, protease inhibitor cocktail

Host cell: HEK293T

Transfection: Expression vector containing full-length human RXFP2 cDNA (GenBank accession number: NM_021634)

Recommended Storage: Store at -80°C. Avoid repeated freeze-thaw.

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.

Functional Assay:

Figure 1

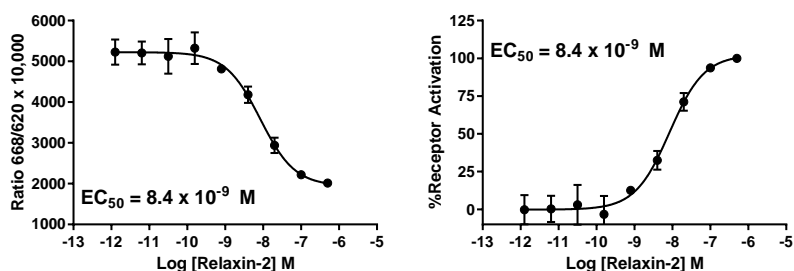


Figure 1. Dose-dependent increase of intracellular cAMP level upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC) using RXFP2 stable cells.

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.

© 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.