

MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT RXFP1 RECEPTOR

PRODUCT INFORMATION

Catalog Number: MC1287NTa

Lot Number: MC1287NTa-010516

Quantity: 1 vial (10.8mg/mL, 1mg)

Packaging Buffer: 20mM Gly-Gly,
1mM MgCl₂, 25 mM Sucrose (pH 7.2)

Host cell: HEK293T

Transfection: Expression vector
containing full-length human RXFP1
cDNA (GenBank accession number:
NM_021634)

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

Data sheet

Background: The relaxin family peptide receptor RXFP1, also known as LGR7 or RX1, is a 757-amino acid 7 transmembrane protein. RT-PCR detected RXFP1 expression in various organs of the reproductive system along with the brain, heart, and lungs. RXFP1 is known to be important to the connective tissue cells of the breast and play a central role in endometrial decidualization.

Application: Functional assays

Figure 1

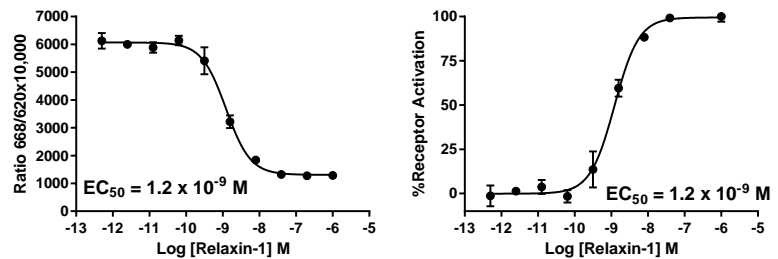


Figure 2

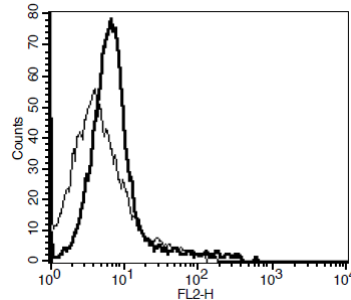


Figure 1. Dose-dependent increase of intracellular cAMP level upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC) using RXFP1 stable cells. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-RXFP1 antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Ivell *et al.* (2003) Immunoexpression of the relaxin receptor LGR7 in breast and uterine tissues of humans and primates. *Reprod Biol Endocrinol* 1:114-114.

Hsu *et al.* (2002) Activation of orphan receptors by the hormone relaxin. *Science* 295:671-674.

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.