

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT RXFP1 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1287

Lot Number: DC1287-073114

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human RXFP1 cDNA (GenBank accession number: NM_021634) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS

Stability: Stable for 1-2 days after thawing

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

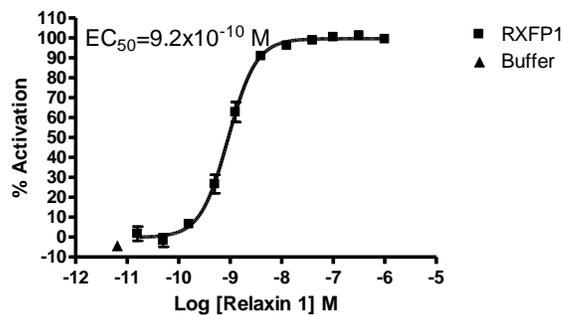


Figure 1. Dose-dependent increase of intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

References:

Ivell *et al.* (2003) Immunoeexpression 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.

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.