

**MULTISCREEN™ STABLE CELL LINE**  
**HUMAN RECOMBINANT RXFP1 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** C1287NTb

**Lot Number:** C1287NTb-091015

**Quantity:** 1 vial ( $2 \times 10^6$ ) 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)

**Recommended Storage:** Liquid nitrogen upon receiving

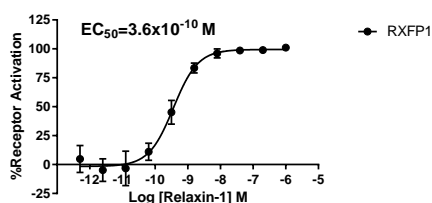
**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ g/mL puromycin

**Stability:** Stable after minimum of two months continuous growth

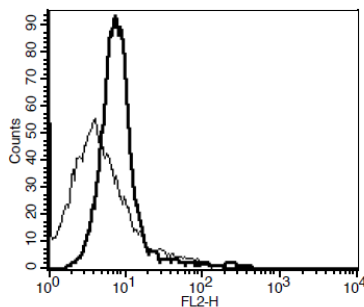
**Background:** The relaxin family peptide receptor RXFP1, also known as LGR7 or RX1, is a 757-amino acid 7 transmembrane protein. RT-PCR detected RXFP1 expression in various organs of the reproductive system along with the brain, heart, and lungs. RXFP1 is known to be important to the connective tissue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

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