

## MULTISCREEN™ STABLE CELL LINE MEMBRANE RECOMBINANT HUMAN RXFP1 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** MC1287

**Lot Number:** MC1287-04162013

**Quantity:** 1 vial (14.78 mg/mL, 5 mg)

**Packaging Buffer:** 20 mM Gly-Gly, 1 mM MgCl<sub>2</sub>, 25 mM Sucrose (pH 7.2)

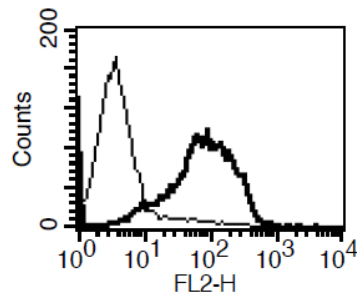
**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human RXFP1 cDNA (GenBank accession number: NM\_021634) with FLAG tag sequence at N-terminus

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

**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.

**Figure 1**



**Figure 1.** Histogram showing receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG 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.