

## MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT RXFP3 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** MC1306

**Lot Number:** MC1306-032414

**Quantity:** 13.9 mg/ml, 10 mg

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

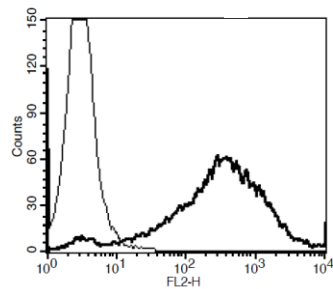
**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human RXFP3 cDNA (GenBank accession number: NM\_016568.2) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

**Background:** The relaxin family peptide receptor RXFP3, also known as GPCR135, RLN3R1 or SALPR, is a 469-amino acid, 7-transmembrane protein. RT-PCR detected RXFP3 expression in various sections of the brain and the testis. RXFP3 is known to be important to several functions of the central nervous system, especially sensory processing under stressful conditions.

**Figure 1**



**Figure 1.** 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:

Cgen *et al.* (2005) Pharmacological characterization of relaxin-3/INSL7 receptors GPCR135 and GPCR142 from different mammalian species. *J Pharmacol Exp Ther* 312:83-95.

Liu *et al.* (2003) Identification of relaxin-3/INSL7 as an endogenous ligand for orphan G-protein-coupled receptor GPCR135. *J Biol Chem* 278:50754-50764.

**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.