

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT CXCR7 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1150

**Lot Number:** C1150-040809

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

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** HEK293T

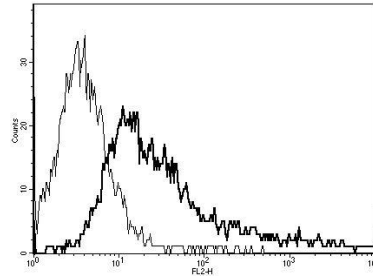
**Transfection:** Expression vector containing full-length human CXCR7 cDNA (GenBank Accession Number: NM\_020311.2) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

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

**Stability:** in progress

**Background:** CXCR7 (or RDC1) is a recently orphanized G-protein coupled receptor which binds with high affinity the inflammatory and homing chemokines CXCL11/ITAC and CXCL12/SDF-1. CXCR7 is expressed in bladder, spleen, heart, skeletal muscle, peripheral nervous system and placenta. CXCR7 does not mediate typical chemokine receptor responses such as leukocyte trafficking. Recent findings in zebrafish indicate that a critical activity of the receptor is scavenging of CXCL12 thereby generating guidance cues for CXCR4-dependent migration.



**Figure Legend:** 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:

Sreedharan *et al.* (1991) Cloning and expression of the human vasoactive intestinal peptide receptor. *Proc Natl Acad Sci USA* 88:4986-4990.

Shimizu *et al.* (1991) A putative G protein-coupled receptor, RDC1, is a novel coreceptor for human and simian immunodeficiency viruses. *Virology* 74:619-626.

**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.