

MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT CXCR7 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1150

Lot Number: C1150-040809

Quantity: 1 vial (2×10^6) 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 μ 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.

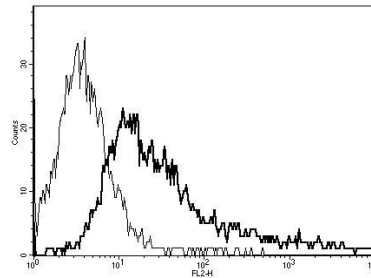


Figure Legend: Receptor Expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

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.

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