

MULTISCREEN™ STABLE CELL LINE MEMBRANE RECOMBINANT HUMAN CHEMOKINE CXCR4 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: MC1004

Lot Number: 07/15/11

Quantity: 1 vial (11.8 mg/mL, 1 mg)

Packaging Buffer: 20 mM Gly-Gly, 1 mM MgCl₂, protease 25 mM Sucrose (pH 7.2)

Host cell: HEK293T

Transfection: Expression vector containing full-length human CXCR4 cDNA (GenBank Accession Number NM_003467) with FLAG tag sequence at N-terminus

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

Background: CXCR4 is a receptor for the C-X-C chemokine SDF-1 (Stromal Cell-Derived Factor 1). It is involved in haematopoiesis and cardiac ventricular septum formation, and plays an essential role in vascularization of the gastrointestinal tract, cerebellar development and survival of hippocampal-neuron of central nerve system. CXCR4 also acts as a primary receptor for some HIV-2 isolates and as a co-receptor with CD4 for HIV-1 X4 viruses.

Figure 1

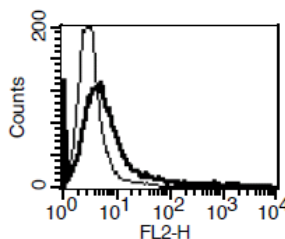


Figure 1. Histogram showing receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Hernandez *et al.* (2003) Mutations in the chemokine receptor gene CXCR4 are associated with WHIM syndrome, a combined immunodeficiency disease. *Nat Genet* 34:70-74.

Juarez *et al.* (2004) Chemokines and their receptors as therapeutic targets: the role of the SDF-1/CXCR4 axis. *Curr Pharm Des* 10:1245-1259.

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.