

## MULTISCREEN™ STABLE CELL LINE MEMBRANE RECOMBINANT HUMAN CHEMOKINE CXCR4 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** MC1004-1

**Lot Number:** 07/15/11

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

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

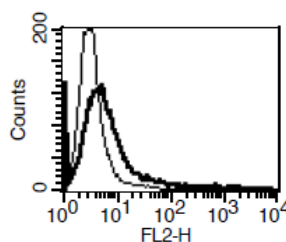
**Host cell:** HEK293T

**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.**

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