

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT XCR1 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1007

Lot Number: C1007-042710

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 XCR1 cDNA (GenBank accession number NM_005283.2) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μ g/mL puromycin

Stability: Stable in culture for minimum of two months

Background: Chemokines and their receptors play a vital role in the regulation of responses in the immune system and in the recruitment of certain lymphocytes. The human chemokine receptor XCR1 is a G-protein-coupled receptor with 333 amino acids. Specifically, XCR1 has been associated with the induction of chemotaxis and has been identified in neutrophils, t-lymphocytes, b-lymphocytes through RT-PCR.

Application: Functional assays

Figure 1

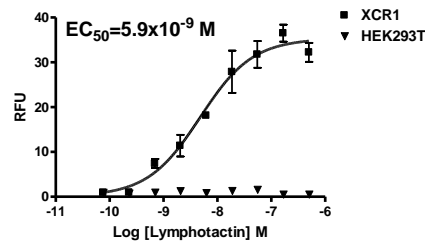


Figure 2

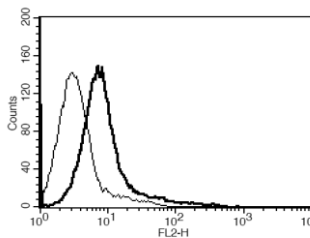


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Huang *et al.* (2001) Neutrophils and B cells express XCR1 receptor and chemotactically respond to lymphotoctin. *Biochem Biophys Res Commun* 281:378-382.

Yoshida *et al.* (1998) Identification of single C motif-1/lymphotoctin receptor XCR1. *J Biol Chem* 273:16551-16554.

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.