

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT XCR1 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** C1007

**Lot Number:** C1007-042710

**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 XCR1 cDNA (GenBank accession number NM\_005283.2) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

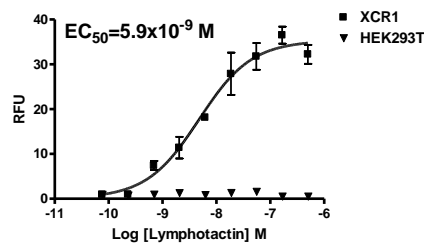
**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ g/mL puromycin

**Stability:** Stable in culture for minimum of two months

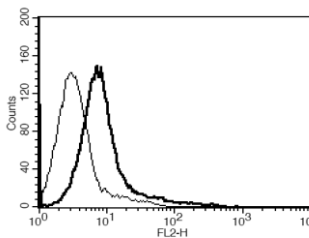
**Background:** Chemokines and their receptors play a vital role in the regulation of responses in the immune system and in the recruitment of certain lymphocytes. The human chemokine receptor XCR1 is a G-protein-coupled receptor with 333 amino acids. Specifically, XCR1 has been associated with the induction 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, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **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 lymphotactin. *Biochem Biophys Res Commun* 281:378-382.

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

**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.