

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

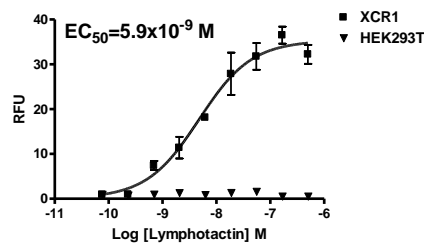


Figure 2

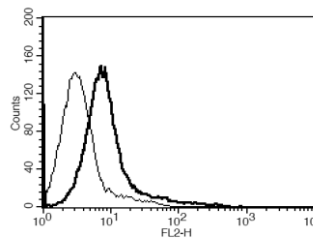


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

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.