

MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT CXCR3 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: CG1003

Lot Number: CG1003-030612

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T Gαq5

Transfection: Expression vector containing full-length human CXCR3 cDNA (GenBank Accession Number: NM_001504) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μg/mL puromycin, 250 μg/ml hygromycin

Stability: In progress

Background: CXCR3 (chemokine receptor 3) is a receptor for SCYB9/MIG, SCYB10/INP10 and SCYB11/ITAC. The chemokines that target CXCR3 in concert with Th1 cytokine appear to play a beneficial role to atopic disorders including a range of conditions such as allergic asthma, -rhinitis, -conjunctivitis, -dermatitis, food and drug allergies and anaphylaxis.

Application: Functional assays

Figure 1

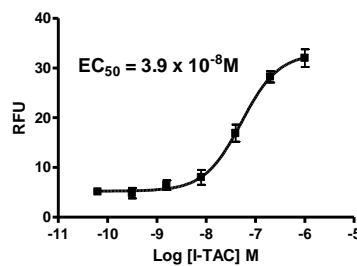


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Loetscher, M., B. Gerber, et al. (1996). "Chemokine receptor specific for IP10 and mig: structure, function, and expression in activated T-lymphocytes." *J Exp Med* 184(3): 963-9.

Arimilli, S., W. Ferlin, et al. (2000). "Chemokines in autoimmune diseases." *Immunol Rev* 177: 43-51.

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.