

**MULTISCREEN™ STABLE CELL LINE**  
**HUMAN RECOMBINANT CXCR3 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** CG1003

**Lot Number:** CG1003-030612

**Quantity:** 1 vial ( $2 \times 10^6$ ) 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

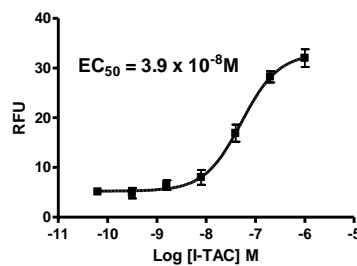
**Propagation Medium:** DMEM, 10% FBS, 1 µg/mL puromycin, 250 µg/ml hygromycin

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

**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.