

## MULTISCREEN™ MEMBRANE PREPARATION HUMAN RECOMBINANT CXCR3 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** MCG1003

**Lot Number:** MCG1003-050614

**Quantity:** 11.8 mg/ml, 1 mg

**Packaging Buffer:** 20 mM Gly-Gly, 1 mM MgCl<sub>2</sub>, 25 mM Sucrose (pH 7.2)

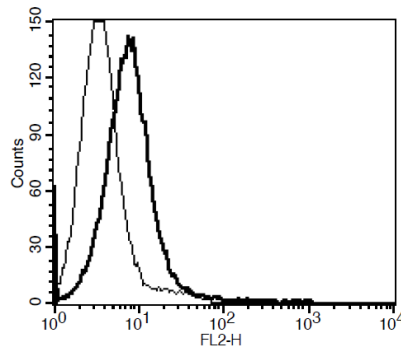
**Host cell:** HEK293T Gaqi5

**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

**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.

**Figure 1**



**Figure 1.** 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:

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.**

© 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.