

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT CXCR5 RECEPTOR

### PRODUCT INFORMATION

**Catalog Number:** C1005-1a

**Lot Number:** C1005-1a-062210

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

**Freeze Medium:** Sigma Freezing Medium (C-6164)

**Host cell:** CHO-K1

**Transfection:** Expression vector containing full-length human CXCR5 cDNA (GenBank Accession Number NM\_001716.2) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM/F12, 10% FBS, 10  $\mu$ g/mL puromycin

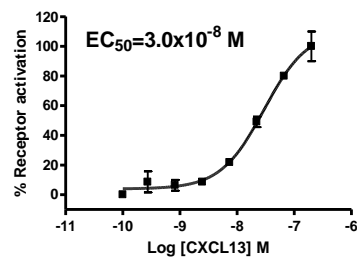
**Stability:** Stable in culture for a minimum of 1 month

### Data sheet

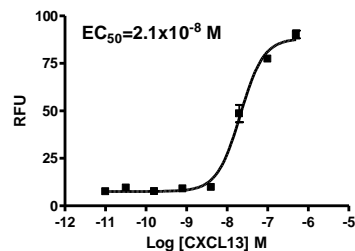
**Background:** CXCR5 (CXC-chemokine receptor 5, or Burkitt lymphoma receptor 1) is a receptor for CXCL13 and expresses in mature B-cells, a subpopulation of T-cells and Burkitt lymphoma cells. CXCR5 exerts possibly a regulatory function in Burkitt lymphoma lymphomagenesis and/or B-cell differentiation. It is a potential candidate for cell-cell interaction and activation of mature B-lymphocytes in lymphatic tissues.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC).

**Figure 2.** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation.

### References:

Dobner *et al.* (1992) Differentiation-specific expression of a novel G protein coupled receptor from Burkitt's lymphoma. *Eur J Immunol* 22:2795-2799.

Campbell *et al.* (2003) Chemokines in the systemic organization of immunity. *Immunol Rev* 195:58-71.

**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.