

**MULTISCREEN™ STABLE CELL LINE  
HUMAN RECOMBINANT CCR5 RECEPTOR**

**PRODUCT INFORMATION**

**Catalog Number:** C1013a

**Lot Number:** C1013a-050710

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

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

**Host cell:** HEK293T

**Transfection:** Full-length human CCR5 cDNA (GenBank Accession Number NM\_000579) with flag tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ g/mL puromycin

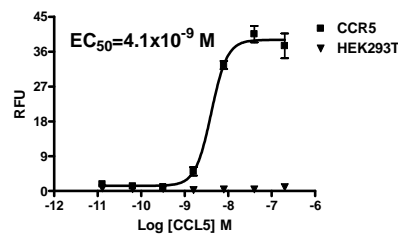
**Stability:** Stable in culture for minimum of two months

**Data sheet**

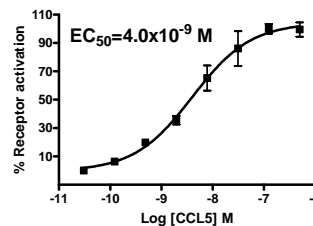
**Background:** CCR5 (chemokine (C-C motif) receptor 5) is a receptor for C-C type chemokines MIP-1- $\alpha$ , MIP-1- $\beta$  and CCL5. CCR5 plays a role in the control of granulocytic lineage proliferation or differentiation. It also acts as co-receptor with CD4 for primary non-syncytium-inducing strains (macrophage-tropic) of HIV-1 virus.

**Application:** Functional assays

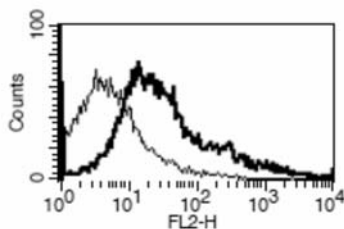
**Figure 1**



**Figure 2**



**Figure 3**



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP accumulation upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC). **Fig 3.** 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:**

Deng *et al.* (1996) Identification of a major co-receptor for primary isolates of HIV-1. *Nature* 381:661-666.

Oppermann (2004) Chemokine receptor CCR5: insights into structure, function, and regulation. *Cell Signal* 16:1201-1210.

**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.