

**MULTISCREEN™ DIVISION ARRESTED CELL LINE  
HUMAN RECOMBINANT CCR5 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1013b

**Lot Number:** 03/12/13

**Quantity:** 1 vial ( $4 \times 10^6$ ) frozen cells

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

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human CCR5 cDNA (GenBank accession number NM\_000579) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

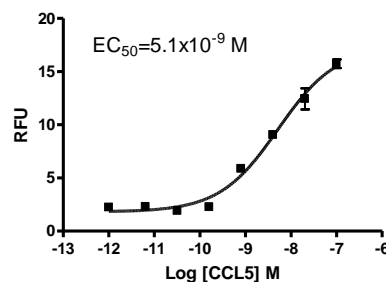
**Propagation Medium:** DMEM, 10% FBS

**Stability:** Stable for 1-2 days after thawing

**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**



Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). Cells were transiently transfected with Gqj5.

**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.**

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.