

MULTISCREEN™ DIVISION-ARRESTED CELL LINE
HUMAN RECOMBINANT CCR5 RECEPTOR

PRODUCT INFORMATION

Catalog Number: DC1013a

Lot Number: DC1013a-032417

Quantity: 1 vial (4×10^6) 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

Stability: Stable for 1-2 days after thawing

Data sheet

Background: CCR5 (chemokine (C-C motif) receptor 5) is a receptor for C-C type chemokines MIP-1- α , MIP-1- β 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

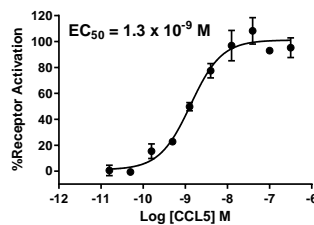


Figure 1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP accumulation upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

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.

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