

**MULTISCREEN™ DIVISION ARRESTED CELL LINE  
HUMAN RECOMBINANT CCR4 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1012

**Lot Number:** 7/19/11

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

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

**Host cell:** HEK293T

**Transfection:** Expression vector containing full-length human CCR4 cDNA (GenBank accession number NM\_005508.2) with FLAG tag sequence at N-terminus

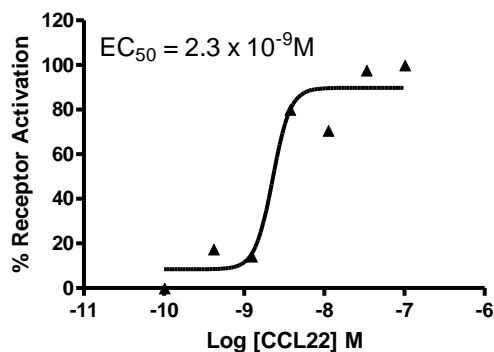
**Recommended Storage:** Liquid nitrogen upon receiving

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

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

**Background:** CCR4 (C-C chemokine receptor type 4) is a high affinity receptor for the C-C type chemokines CCL17/TARC and CCL22/MDC. CCR4 receptors express in the thymus, peripheral blood leukocytes and the spleen. CCR4 functions as a chemoattractant homing receptor on circulating memory lymphocytes and as a coreceptor for some primary HIV-2 isolates. In the CNS, CCR4 mediates hippocampal-neuron survival. CCR4 is one of the type 2 T helper cell (Th2) markers. Th2 cells are recruited to the asthmatic lung in response to allergen challenge and CCR4 has been therefore considered as a therapeutic target of lung inflammation, including eosinophilic inflammation of the lung and asthma. In addition, CCR4 has been reported as a potential independent and significant prognostic factor for adult T-cell leukemia/lymphoma (ATLL).

**Application:** cAMP assays



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

**References:**

Power and Meyer (1995) Molecular cloning and functional expression of a novel CC chemokine receptor cDNA from a human basophilic cell line. *J Biol Chem* 270:19495-19500.

Imai and Baba (1997) The T cell-directed CC chemokine TARC is a highly specific biological ligand for CC chemokine receptor 4. *J Biol Chem* 272:15036-15042.

**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.