

**MULTISCREEN™ DIVISION ARRESTED CELL LINE  
HUMAN RECOMBINANT CCR2 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1010

**Lot Number:** 01/16/13

**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 CCR2 cDNA (GenBank accession number NM\_000647.3) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

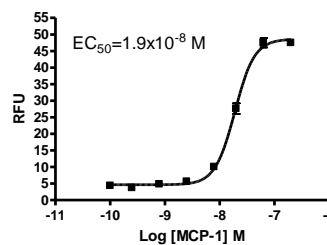
**Propagation Medium:** DMEM, 10% FBS

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

**Background:** CCR2 (C-C chemokine receptor type 2 or monocyte chemoattractant protein 1 receptor) is a receptor for the MCP-1, MCP-3 and MCP-4 chemokines. It is also an alternative coreceptor with CD4 for HIV-1 infection. CCR2 has been widely considered as a potential therapeutic target for autoimmune disease and inflammatory disorders of the lung, particularly rheumatoid arthritis, and various CCR2 blocking agents have been developed, some of which might have therapeutic values.

**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 Gαq15 before division arrest.

**References:**

Quinones *et al.* (2005) The complex role of the chemokine receptor CCR2 in collagen-induced arthritis: implications for therapeutic targeting of CCR2 in rheumatoid arthritis. *J Mol Med* 83:672-681.

Rose *et al.* (2003) Significant involvement of CCL2 (MCP-1) in inflammatory disorders of the lung. *Microcirculation* 10:273-288.

**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.