

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

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1010

Lot Number: 01/16/13

Quantity: 1 vial (4×10^6) 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

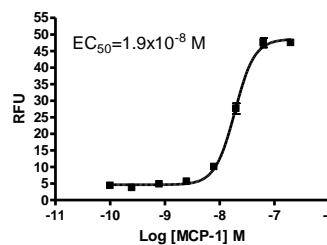
Propagation Medium: DMEM, 10% FBS

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 Gaq15 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.