

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT CCR1 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** CG1009

**Lot Number:** CG1009-02062013

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

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

**Host cell:** HEK293T Gqi5

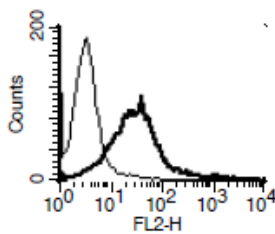
**Transfection:** Expression vector containing full-length human CCR1 cDNA (GenBank Accession Number NM\_001295) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS, 250 µg/mL hygromycin, 1µg/mL puromycin

**Stability:** Stable in culture for minimum of two months

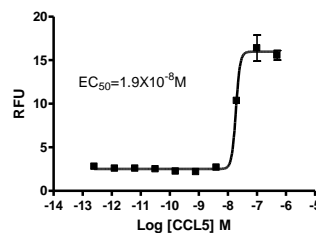
#### Figure 2



**Background:** CCR1 is a CC chemokine receptor with high affinity for RANTES. It is demonstrated that targeted disruption of the CCR1 receptor is associated with protection from pulmonary inflammation secondary to acute pancreatitis in the mouse. The protection from lung injury is associated with decreased levels of TNF-alpha in a temporal sequence indicating that the activation of the CCR1 receptor is an early event in the systemic inflammatory response syndrome.

**Application:** Functional assays

#### Figure 1



**Figure 1.** Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 3.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

#### References:

- Wieser *et al.* (2005) Expression and regulation of CCR1 in peritoneal macrophages from women with and without endometriosis. *Fertil Steril* 83:1878-1881.
- Neote *et al.* (1993) Molecular cloning, functional expression, and signaling characteristics of a C-C chemokine receptor. *Cell* 72:415-425.

**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.