

## MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT CCR8 RECEPTOR

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** CG1016

**Lot Number:** CG1016-042513

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

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

**Host cell:** HEK293T Gq*α*5

**Transfection:** Expression vector containing full-length human CCR8 cDNA (GenBank Accession Number NM\_005201.2) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

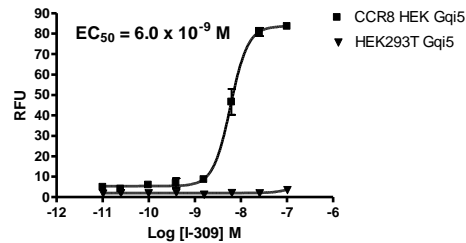
**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ g/mL puromycin, 250  $\mu$ g/mL hygromycin

**Stability:** In progress

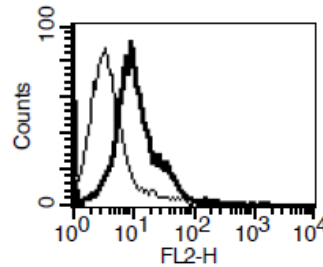
**Background:** CCR8 (C-C chemokine receptor type 8) is a receptor for the chemokines CCL1 (I-309). The receptor regulates monocyte chemotaxis and thymic cell line apoptosis, and is an additional alternative coreceptor with CD4 for HIV-1 infection. CCR8 plays a role on cell types of relevance to respiratory diseases such as asthma, chronic obstructive pulmonary disease (COPD) and chronic bronchitis. The interference of CCR8 from its ligand represents a new approach in allergy therapy.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**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 2.** 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:

Simmons *et al.* (2000) Co-receptor use by HIV and inhibition of HIV infection by chemokine receptor ligands. *Immunol Rev* 177:112-126.

Garcia *et al.* (2005) New chemokine targets for asthma therapy. *Curr Allergy Asthma Rep* 5:155-160

**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.