

**MULTISCREEN™ STABLE CELL LINE  
HUMAN RECOMBINANT CCR7 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** CG1015

**Lot Number:** CG1015-03052013

**Quantity:** 1 vial (2 x 10<sup>6</sup>) frozen cells

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

**Host cell:** HEK293T Gqi5

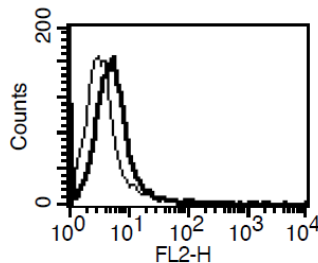
**Transfection:** Expression vector containing full-length human CCR7 cDNA (GenBank Accession Number: NM\_001838) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

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

**Stability:** In progress

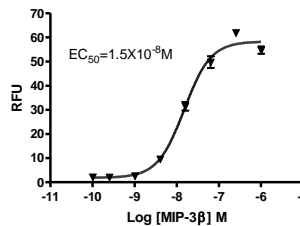
**Figure 2**



**Background:** CCR7 (C-C chemokine receptor type 7) is a receptor for the chemokine MIP-3β. CCR7 has been identified as a key regulator of homeostatic B and T cell trafficking to secondary lymphoid organs. CCR7 is also an essential mediator for entry of both dermal and epidermal dendritic cells into the lymphatic vessels within the dermis. CCR7 plays an important role in tumor cell migration and lymph node metastasis. Expression of CCR7 is related to the development of lymph node metastasis in nonsmall cell lung cancers.

**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 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:**

Takanami (2003) Overexpression of CCR7 mRNA in nonsmall cell lung cancer: correlation with lymph node metastasis. *Int J Cancer* 105:186-189.

Ohl *et al.* (2004) CCR7 governs skin dendritic cell migration under inflammatory and steady-state conditions. *Immunity* 21:279-288.

**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.