

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

Data sheet

PRODUCT INFORMATION

Catalog Number: C1015

Lot Number: C1015-023006

Quantity: 1 vial (2×10^6) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

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 μ g/mL puromycin

Stability: Stable in culture for minimum of two months

Background: CCR7 (C-C chemokine receptor type 7) is a receptor for the chemokine MIP-3 β . 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

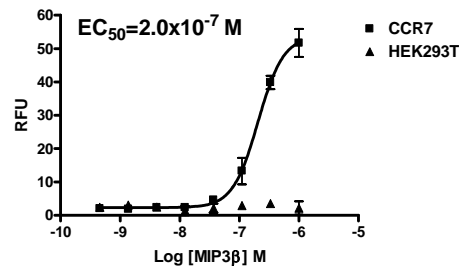


Figure 2

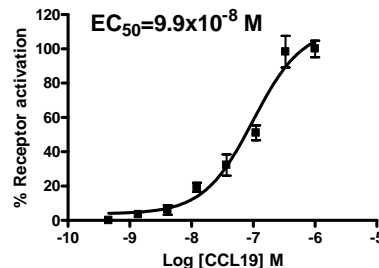


Figure 3

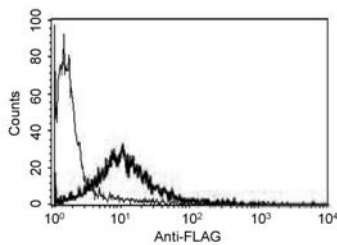


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP accumulation upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC). **Figure 3.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

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.

© 2005 Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.

www.multispaninc.com
sales@multispaninc.com
support@multispaninc.com

Ver. October 2005

Phone: +1 (510) 887-0817
Fax: +1 (510) 887-0863
26219 Eden Landing Road
Hayward, CA 94545-3718
U.S.A.