

**MULTISCREEN™ DIVISION ARRESTED CELL LINE  
HUMAN RECOMBINANT CCR7 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** DC1015

**Lot Number:** 01/16/13

**Quantity:** 1 vial ( $4 \times 10^6$ ) 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

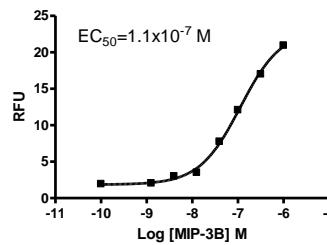
**Propagation Medium:** DMEM, 10% FBS

**Stability:** Stable for 1-2 days after thawing

**Background:** CCR7 (C-C chemokine receptor type 7 is a receptor for the chemokine MIP-3 $\beta$ . 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**



Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. Cells were transiently transfected with G $\alpha$ q15 before division arrest.

**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.