

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

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1015

Lot Number: 01/16/13

Quantity: 1 vial (4×10^6) 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

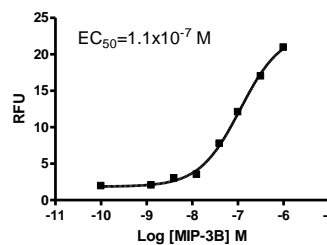
Propagation Medium: DMEM, 10% FBS

Stability: Stable for 1-2 days after thawing

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



Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). Cells were transiently transfected with Gaq15 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.

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