

## MULTISCREEN™ STABLE CELL LINE DOG RECOMBINANT CCR9 RECEPTOR

### PRODUCT INFORMATION

**Catalog Number:** CGd1017

**Lot Number:** CGd1017-111213

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

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

**Host cell:** HEK293T Gαq5

**Transfection:** Expression vector containing full-length human CCR9 cDNA (GenBank Accession Number: NM\_001284476) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

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

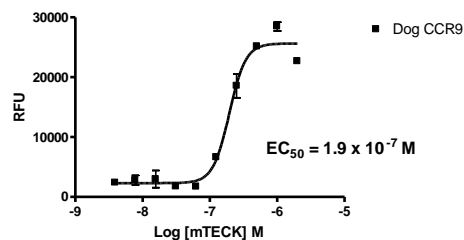
**Stability:** Stable in culture for minimum of two months

### Data sheet

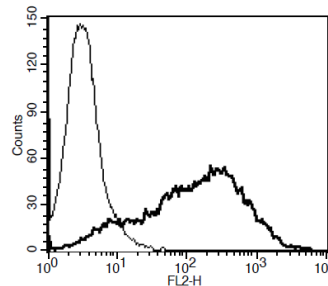
**Background:** CCR9 (C-C chemokine receptor type 9) is a thymus-specific receptor for chemokine CCL25 (SCYA25 or TECK). It is an alternative co-receptor with CD4 for HIV-1 infection. Activation of the receptor may lead to potent cFLIP(L)-independent resistance to cycloheximide-induced apoptosis and modest resistance to Fas-mediated apoptosis possibly via activation of multiple signaling components involving Akt and glycogen synthase kinase 3β. Thus, dissection of signaling components involved in the CCR9-mediated antiapoptosis could be a framework for cell survival mechanisms and may provide options for therapeutic interventions.

**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:

Youn *et al.* (2002) Role of the CC chemokine receptor 9/TECK interaction in apoptosis. *Apoptosis* 7:271-276.

Zaballos *et al.* (1999) Cutting edge: identification of the orphan chemokine receptor GPR-9-6 as CCR9, the receptor for the chemokine TECK. *J Immunol* 162(10):5671-5.

**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.