

**MULTISCREEN™ STABLE CELL LINE  
HUMAN RECOMBINANT CCR6 RECEPTOR**

**Data Sheet**

**PRODUCT INFORMATION**

**Catalog Number:** C1014b

**Lot Number:** C1014b-031516

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

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

**Host cell:** HEK293T

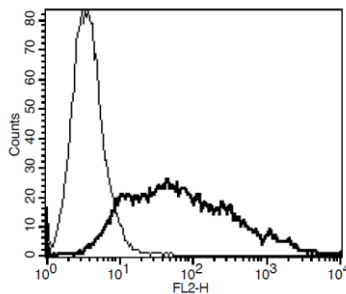
**Transfection:** Expression vector containing full-length human CCR6 cDNA (GenBank Accession Number NM\_004367) with FLAG tag sequence at N-terminus

**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ g/mL puromycin

**Stability:** Stability in progress

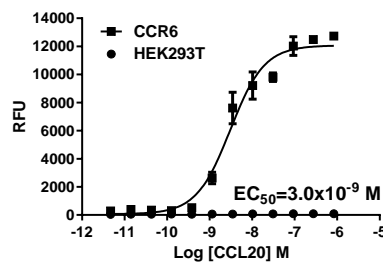
**Figure 3**



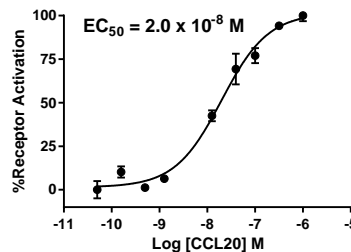
**Background:** CCR6 is a receptor for macrophage inflammatory protein (MIP)-3 $\alpha$  (CCL20). It is expressed in immature dendritic cells as well as in B lymphocytes and memory T cells. In gene-targeted mice lacking CCR6, myeloid dendritic cells are either absent or underrepresented in the subepithelial dome compared with wild-type mice. CCR6 and its ligands contribute to the normal homeostatic regulation of extrathymic T cell development in the intestine.

**Application:** Functional assays

**Figure 1**



**Figure 2**



**Figure 1.** Dose-dependent calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01). **Figure 2.** Dose-dependent inhibition of forskolin-stimulated intracellular cAMP accumulation upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **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:**

Baba et al. (1997) Identification of CCR6, the specific receptor for a novel lymphocytedirected CC chemokine LARC. *J Biol Chem* 272:14893-14898.

Cook et al. (2000) CCR6 mediates dendritic cell localization, lymphocyte homeostasis, and immune responses in mucosal tissue. *Immunity* 12:495-503.

Lugering et al. (2003) Lymphoid precursors in intestinal cryptopatches express CCR6 and undergo dysregulated development in the absence of CCR6. *J Immunol* 171:2208-2215.

**FOR RESEARCH USE ONLY.**