

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT CXCR6 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: C1006-1

Lot Number: C1006-1-062210

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: CHO-K1

Transfection: Full-length Human CXCR6 cDNA (GenBank Accession Number NM_006564) with FLAG-tag sequence at the N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS, 10 μ g/mL puromycin

Stability: Stable in culture for minimum of two months

Background: CXCR6 (CXC-chemokine receptor type 6) is a receptor for CXC chemokine CXCL16. CXCR6 is expressed by naive CD8 T cells, natural killer T cells, a subset of memory CD4 T cells and plasma cells. CXCR6 and its ligand play an important role in the trafficking of the cells expressing them. It is demonstrated that CXCL16⁺ macrophages in synovial fluid of rheumatoid arthritis patients lead to recruitment of CXCR6⁺ memory T cells, thereby contributing to the inflammatory cascade associated with rheumatoid arthritis pathology. CXCR6 is also used as a coreceptor by simian immunodeficiency viruses (SIVs) and by strains of HIV-2 and M-tropic HIV-1.

Application: Functional assays

Figure 1

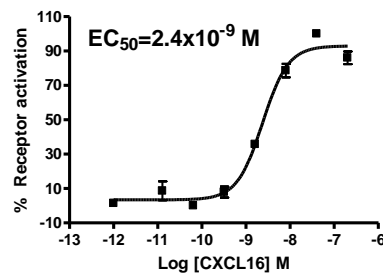


Figure 2

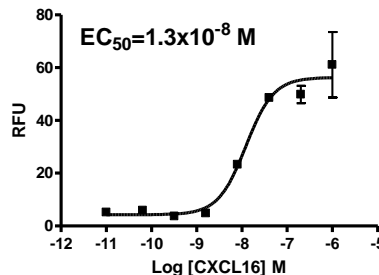


Figure 1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP level upon treatment with ligand, measured with Multiscreen™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01). **Figure 2.** Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™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

References:

Deng *et al.* (1997) Expression cloning of new receptors used by simian and human immunodeficiency viruses. *Nature* 388:296-300.

Blaak *et al.* (2005) CCR5, GPR15, and CXCR6 are major coreceptors of human immunodeficiency virus type 2 variants isolated from individuals with and without plasma viremia. *J Virol* 79:1686-1700.

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.