

MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT CXCR6 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1006

Lot Number: C1006-070711

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

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, 10% FBS, 1 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

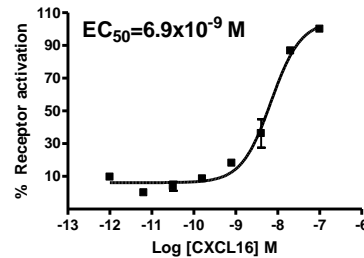
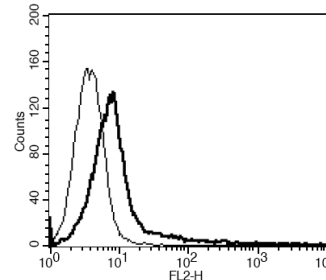


Figure 2



1. Dose-dependent inhibition of forskolin-stimulated intracellular cAMP accumulation upon treatment with ligand, measured with cAMP HiRange kit (Cisbio 62AM6PEC). **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:

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.

© 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.