

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT CXCR5 RECEPTOR

PRODUCT INFORMATION

Catalog Number: C1005

Lot Number: C1005-060909

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Expression vector containing full-length human CXCR5 cDNA (GenBank Accession Number NM_001716.2) with FLAG tag sequence at 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

Data sheet

Background: CXCR5 (CXC-chemokine receptor 5, or Burkitt lymphoma receptor 1) is a receptor for CXCL13 and expresses in mature B-cells, a subpopulation of T-cells and Burkitt lymphoma cells. CXCR5 exerts possibly a regulatory function in Burkitt lymphoma lymphomagenesis and/or B-cell differentiation. It is a potential candidate for cell-cell interaction and activation of mature B-lymphocytes in lymphatic tissues.

Application: Functional assays

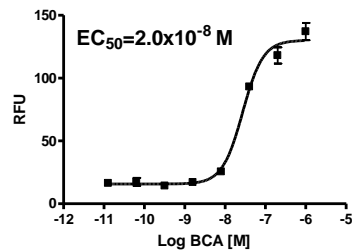


Figure legend: Dose-dependent calcium flux upon treatment with ligand, monitored with FlexStation.

References:

Dobner *et al.* (1992) Differentiation-specific expression of a novel G protein coupled receptor from Burkitt's lymphoma. *Eur J Immunol* 22:2795-2799.

Campbell *et al.* (2003) Chemokines in the systemic organization of immunity. *Immunol Rev* 195:58-71.

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.