

**MULTISCREEN™ DIVISION ARRESTED CELL LINE
HUMAN RECOMBINANT CXCR5 RECEPTOR**

Data sheet

PRODUCT INFORMATION

Catalog Number: DC1005-1a

Lot Number: 011818

Quantity: 1 vial (4×10^6) frozen cells

Freeze Medium: Cellbanker 2 (Amsbio 11891)

Host cell: CHO-K1

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

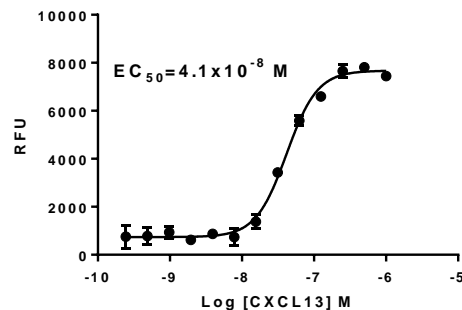
Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM/F12, 10% FBS

Stability: Stable for 1-2 days after thawing

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



Dose-dependent stimulation of calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

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.

Multispan Inc. All rights reserved. No part of this document may be reproduced in any form without prior permission in writing.