

MULTISCREEN™ STABLE CELL LINE HUMAN RECOMBINANT CMKLR1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: CG1147

Lot Number: CG1147-022510

Quantity: 1 vial (2 x 10⁶) frozen cells

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T Gaqi5

Transfection: Expression vector containing full-length human CMKLR1 cDNA (GenBank Accession Number: NM_004072) with FLAG tag sequence at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 µg/mL puromycin, 150 µg/mL hygromycin

Stability: Stable after minimum of two months continuous growth

Background: Chemokine receptor-like 1 (CMKLR1) (or ChemR23) is a receptor for TIG2 and Chemerin. ChemR23/Chemerin may play a key role in directing plasmacytoid dendritic cells trafficking. The receptor also acts as a coreceptor for several SIV strains (SIVMAC316, SIVMAC239, SIVMAC17E-FR and SIVSM62A), as well as a primary HIV-1 strain (92UG024-2).

Application: Functional assays

Figure 1

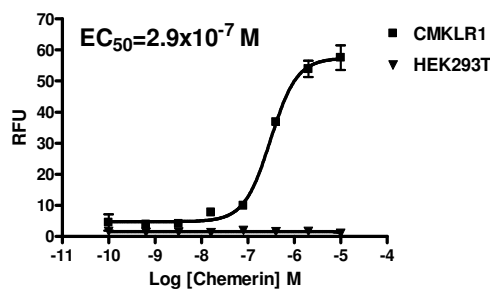


Figure 2

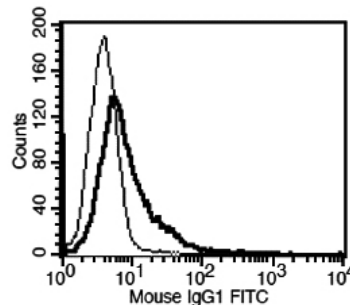


Figure 1. Dose-dependent stimulation of calcium flux upon treatment with ligand, monitored with FlexStation. **Figure 2.** Receptor expression on cell surface measured by flow cytometry (FACS) using an anti-FLAG antibody. Thin line: parental cells; thick line: receptor-expressing cells.

References:

Meder *et al.* (2003) Characterization of human circulating TIG2 as a ligand for the orphan receptor ChemR23. *FEBS Lett* 555:495-499.

Wittamer *et al.* (2004) The C-terminal nonapeptide of mature chemerin activates the chemerin receptor with low nanomolar potency. *J Biol Chem* 279:9956-9962.

Samson *et al.* (1998) ChemR23, a putative chemoattractant receptor, is expressed in monocyte-derived dendritic cells and macrophages and is a coreceptor for SIV and some primary HIV-1 strains. *Eur J Immunol* 28:1689-1700.

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.