

MULTISCREEN™ β -ARRESTIN2 STABLE CELL LINE HUMAN RECOMBINANT CXCR7 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: CA1150-BA2

Lot Number: CA1150-BA2-080921

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

Freeze Medium: Cellbanker 2
(Amsbio)

Host cell: HEK293T β -Arrestin2

Transfection: Expression vector containing full-length human CXCR7 cDNA (GenBank Accession Number: NM_020311.2) with FLAG tag sequence at N-terminus and ARRB2 cDNA (GenBank Accession Number NM_004313.3)

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 1 μ g/mL puromycin, 50 μ g/mL hygromycin

Stability: In progress

Background: CXCR7 (or RDC1) is a recently orphanized G-protein coupled receptor which binds with high affinity the inflammatory and homing chemokines CXCL11/ITAC and CXCL12/SDF-1. CXCR7 is expressed in bladder, spleen, heart, skeletal muscle, peripheral nervous system and placenta. CXCR7 does not mediate typical chemokine receptor responses such as leukocyte trafficking. Recent findings in zebrafish indicate that a critical activity of the receptor is scavenging of CXCL12 thereby generating guidance cues for CXCR4-dependent migration.

Figure 1.

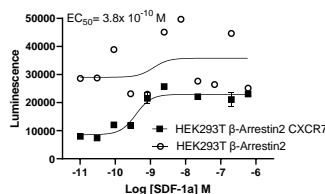


Figure 1. Dose-dependent stimulation from arrestin recruitment upon treatment with ligand, measured with MultiScreen™ β -Arrestin Assay Kit (Multispan MSBA01).

References:

Sreedharan *et al.* (1991) Cloning and expression of the human vasoactive intestinal peptide receptor. *Proc Natl Acad Sci USA* 88:4986-4990.

Shimizu *et al.* (1991) A putative G protein-coupled receptor, RDC1, is a novel coreceptor for human and simian immunodeficiency viruses. *Virology* 74:619-626.

FOR RESEARCH USE ONLY.

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