

MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT CysLT1 RECEPTOR

Data sheet

PRODUCT INFORMATION

Catalog Number: C1308

Lot Number: C1308-073005

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 CysLTR1 cDNA (GenBank Accession Number NM_006639) with FLAG tag at N-terminus

Recommended Storage: Liquid nitrogen upon receiving

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

Stability: Stable after minimum two months continuous growth.

Background: CysLTR1 (Cysteinyl leukotriene receptor 1) is also known as LTD4 receptor (leukotriene D4 receptor), HG55 or HMTMF81. It is a receptor for cysteinyl leukotrienes and has highest affinity to leukotriene D4 (LTD4). The receptor mediates contraction and proliferation of smooth muscle, edema, eosinophil migration and damage to the mucus layer in the lung caused by LTD4. Several antagonists, such as montelukast (Singulair), zafirlukast (Accolate) and pranlukast (Onon), are used in the treatment of asthma.

Application: Functional assays

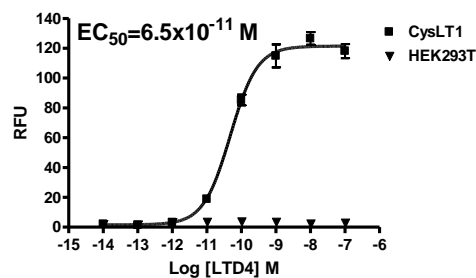


Figure legend: Dose-dependent calcium flux upon treatment with ligand, monitored with FlexStation.

References:

Crowther and Rees (2000) Current treatment of asthma--focus on leukotrienes. *Expert Opin Pharmacother* 1:1021-1040.

Tomari *et al.* (2003) Pranlukast, a cysteinyl leukotriene receptor 1 antagonist, attenuates allergen-specific tumour necrosis factor alpha production and nuclear factor kappa B nuclear translocation in peripheral blood monocytes from atopic asthmatics. *Clin Exp Allergy* 33:795-801.

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.