

**MULTISCREEN™ STABLE CELL LINE**  
**HUMAN RECOMBINANT CysLT1 RECEPTOR**

**Data sheet**

**PRODUCT INFORMATION**

**Catalog Number:** C1308

**Lot Number:** C1308-073005

**Quantity:** 1 vial ( $2 \times 10^6$ ) 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

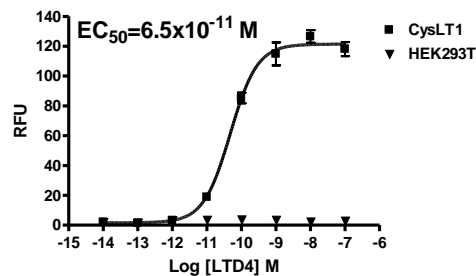
**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS, 1  $\mu$ 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, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

**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.**

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