

**MULTISCREEN™ STABLE CELL LINE
HUMAN RECOMBINANT CysLT2 RECEPTOR**

PRODUCT INFORMATION

Catalog Number: C1307

Lot Number: C1307-082509

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

Freeze Medium: Sigma Freezing Medium (C-6164)

Host cell: HEK293T

Transfection: Full-length human CYSLTR2 cDNA (GenBank Accession Number NM_020377)

Recommended Storage: Liquid nitrogen upon receiving

Propagation Medium: DMEM, 10% FBS, 200 μ g/mL hygromycin

Stability: Stable after minimum two months continuous growth

Data sheet

Background: Cysteinyl leukotriene receptor 2 (CysLTR2) is a receptor for cysteinyl leukotrienes. There are two receptors for cysteinyl leukotrienes, CysLTR1 and CysLTR2. CysLTRs have a clear role in pathophysiological conditions such as asthma, allergic rhinitis and other nasal allergies, and have been implicated in a number of inflammatory conditions in cardiovascular and gastrointestinal diseases.

Application: Functional assays

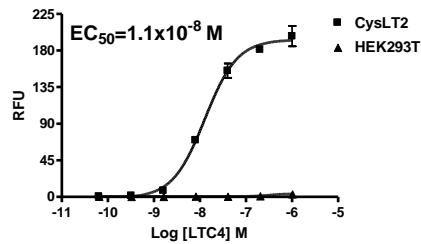


Figure legend: Dose-dependent calcium flux upon treatment with ligand, measured with Multiscreen™ Calcium 1.0 No Wash Assay Kit (Multispan MSCA01).

References:

Takasaki *et al.* (2000) The molecular characterization and tissue distribution of the human cysteinyl leukotriene CysLT(2) receptor. *Biochem Biophys Res Commun* 274:316-322.

Capra (2004) Molecular and functional aspects of human cysteinyl leukotriene receptors. *Pharmacol Res* 50:1-11.

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.