

PRODUCT INFORMATION



CysLT₁ Receptor Polyclonal Antibody Item No. 120500

Overview and Properties

Contents:	This vial contains 500 µl peptide affinity-purified polyclonal antibody.
Synonym:	Cysteinyl Leukotriene Receptor 1
Immunogen:	Synthetic peptide from the C-terminal region of human CysLT ₁
Species Reactivity:	(+) Human, murine; other species not tested
Uniprot No.:	Q9Y271
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Application:	Western blot; the recommended starting dilution 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

The cysteinyl leukotrienes (CysLTs; LTC₄, LTD₄, and LTE₄) contract airway and pulmonary vascular smooth muscle, increase vascular permeability, and stimulate mucus secretion, thereby playing a major role in asthma.¹⁻⁴ LTC₄, LTD₄, and LTE₄ mediate their actions *via* at least two receptors designated CysLT₁ and CysLT₂.¹ Cloning of the human CysLT₁ receptor reveals it is a 337 amino acid protein with a calculated molecular mass of 38,549.⁵ The rank order of binding for leukotrienes to the cloned receptor, as determined using a radioligand binding assay, is LTD₄ >> LTC₄ = LTE₄ >> LTB₄.⁵ The mRNA for the human CysLT₁ receptor is expressed in spleen and peripheral blood leukocytes with smaller amounts in lung, placenta, and small intestine. Immunohistochemical studies reveal the receptor is expressed on eosinophils, mast cells, macrophages, neutrophils, and vascular endothelial cells in the nasal mucosa.⁶

References

1. Gorenne, I., Norel, X., and Brink, C. Cysteinyl leukotriene receptors in the human lung: What's new? *Trends Pharmacol. Sci.* **17(10)**, 342-343 (1996).
2. Dahlén, S.-E., Hansson, G., Hedqvist, P., *et al.* Allergen challenge of lung tissue from asthmatics elicits bronchial contraction that correlates with the release of leukotrienes C₄, D₄, and E₄. *Proc. Natl. Acad. Sci. USA* **80(6)**, 1712-1716 (1983).
3. Busse, W.W. The role of leukotrienes in asthma and allergic rhinitis. *Clin. Exp. Allergy* **26(8)**, 868-879 (1996).
4. Hedqvist, P., Dahlén, S.-E., Gustafsson, L., *et al.* Biological profile of leukotrienes C₄ and D₄. *Acta Physiol. Scand.* **110(3)**, 331-333 (1980).
5. Lynch, K.R., O'Neill, G.P., Liu, Q., *et al.* Characterization of the human cysteinyl leukotriene CysLT₁ receptor. *Nature* **399(6738)**, 789-793 (1999).
6. Shirasaki, H., Kanaizumi, E., Watanabe, K., *et al.* Expression and localization of the cysteinyl leukotriene 1 receptor in human nasal mucosa. *Clin. Exp. Allergy* **32(7)**, 1007-1012 (2002).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/20/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM