

# PRODUCT INFORMATION



## CysLT<sub>1</sub> Receptor Polyclonal FITC Antibody Item No. 14713

### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl peptide affinity-purified, fluorescein-labeled polyclonal antibody.
<b>Synonym:</b>	Cysteinyl Leukotriene Receptor 1
<b>Immunogen:</b>	Synthetic peptide from the C-terminal region of CysLT <sub>1</sub> receptor
<b>Species Reactivity:</b>	(+) Human and mouse; other species not tested
<b>Uniprot No.:</b>	Q9Y271
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥2 years
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Application:</b>	Flow cytometry; the recommended starting dilution is 20 µl/test. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Description

The cysteinyl leukotrienes (CysLTs; LTC<sub>4</sub>, LTD<sub>4</sub>, and LTE<sub>4</sub>) contract airway and pulmonary vascular smooth muscle, increase vascular permeability, and stimulate mucus secretion, thereby playing a major role in asthma.<sup>2-5</sup> LTC<sub>4</sub>, LTD<sub>4</sub>, and LTE<sub>4</sub> mediate their actions *via* at least two receptors designated CysLT<sub>1</sub> and CysLT<sub>2</sub>.<sup>2</sup> Cloning of the human CysLT<sub>1</sub> receptor reveals it is a 337 amino acid protein with a calculated molecular mass of 38,549.<sup>1</sup> The rank order of binding for leukotrienes to the cloned receptor, as determined using a radioligand binding assay, is LTD<sub>4</sub> >> LTC<sub>4</sub> = LTE<sub>4</sub> >> LTB<sub>4</sub>.<sup>1</sup> The mRNA for the human CysLT<sub>1</sub> 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.<sup>6</sup>

### References

1. Lynch, K.R., O'Neill, G.P., Liu, Q., *et al.* Characterization of the human cysteinyl leukotriene CysLT<sub>1</sub> receptor. *Nature* **399**, 789-793 (1999).
2. Gorenne, I., Norel, X., and Brink, C. Cysteinyl leukotriene receptors in the human lung: What's new? *Trends Pharmacol. Sci.* **17**, 342-343 (1996).
3. 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 C<sub>4</sub>, D<sub>4</sub>, and E<sub>4</sub>. *Proc. Natl. Acad. Sci. USA* **80**, 1712-1716 (1983).
4. Busse, W.W. The role of leukotrienes in asthma and allergic rhinitis. *Clin. Exp. Allergy* **26**, 868-879 (1996).
5. Hedqvist, P., Dahlén, S.-E., Gustafsson, L., *et al.* Biological profile of leukotrienes C<sub>4</sub> and D<sub>4</sub>. *Acta Physiol. Scand.* **110**, 331-333 (1980).
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**, 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, 02/02/2026

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