

PRODUCT INFORMATION

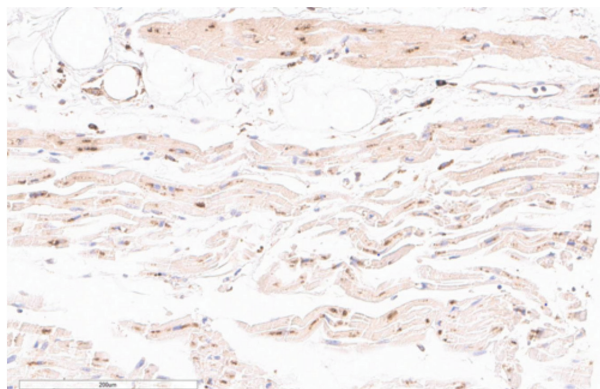


CysLT₂ Receptor (N-Term) Polyclonal Antibody Item No. 120560

Overview and Properties

Contents:	This vial contains peptide affinity-purified polyclonal antibody.
Synonym:	Cysteinyl-Leukotriene Receptor 2 (N-Term)
Immunogen:	Synthetic peptide from the N-terminal region of human CysLT ₂
Cross Reactivity:	(-) CysLT ₁
Species Reactivity:	(+) Human; (-) Rat
Uniprot No.:	Q9NS75
Form:	Lyophilized
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	TBS, pH 7.4, when reconstituted in 500 µl deionized water
Host:	Rabbit
Applications:	ELISA, flow cytometry (FC), immunocytochemistry (ICC), immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution for ELISA is 1:200, ~1:1,000 for FC, and 1:250 for WB. ICC, IHC, and other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human heart tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with CysLT₂ Receptor (N-Term) Polyclonal Antibody (Item No. 120560) at a 1:120 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).

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, 008/14/2019

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

PRODUCT INFORMATION



Description

The cysteinyl leukotrienes (cysLTs) (LTC_4 , LTD_4 , and LTE_4) contract airway and pulmonary vascular smooth muscle, increase vascular permeability and stimulate mucus secretion, thereby playing a major role in asthma.¹⁻⁴ LTC_4 , LTD_4 , and LTE_4 mediate their actions *via* at least 2 receptors designated $CysLT_1$ and $CysLT_2$.¹ Cloning of the human $CysLT_2$ receptor reveals it is a 346 amino acid protein with 38% homology to the human $CysLT_1$ receptor.^{5,6} The rank order of binding for leukotrienes to the cloned receptor, as determined using a radioligand binding assay, is $LTC_4 = LTD_4 \gg LTE_4$.⁵ The mRNA for the human $CysLT_2$ receptor is expressed in lung macrophages, airway smooth muscle, cardiac Purkinje cells, adrenal medulla cells, peripheral blood leukocytes, spleen, placenta, and brain.^{5,7}

References

1. Gorenne, I., Norel, X., and Brink, C. Cysteinyl leukotriene receptors in the human lung: What's new? *Trends Pharmacol. Sci.* **17**, 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 C_4 , D_4 , and E_4 . *Proc. Natl. Acad. Sci. USA* **80**, 1712-1716 (1983).
3. Busse, W.W. The role of leukotrienes in asthma and allergic rhinitis. *Clin. Exp. Allergy* **26**, 868-879 (1996).
4. Hedqvist, P., Dahlén, S.-E., Gustafsson, L., *et al.* Biological profile of leukotrienes C_4 and D_4 . *Acta Physiol. Scand.* **110**, 331-333 (1980).
5. Heise, C.E., O'Dowd, B.F., Figueroa, D.J., *et al.* Characterization of the human cysteinyl leukotriene 2 receptor. *J. Biol. Chem.* **275**, 30531-30536 (2000).
6. Lynch, K.R., O'Neill, G.P., Liu, Q., *et al.* Characterization of the human cysteinyl leukotriene $CysLT_1$ receptor. *Nature* **399**, 789-793 (1999).
7. Takasaki, J., Kamohara, M., Matsumoto, M., *et al.* The molecular characterization and tissue distribution of the human cysteinyl leukotriene $CysLT_2$ receptor. *Biochem. Biophys. Res. Commun.* **274**, 316-322 (2000).

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