

PRODUCT INFORMATION

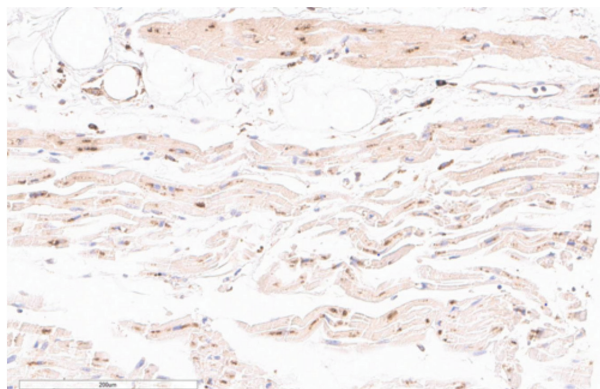


CysLT₂ Receptor (N-Term) Polyclonal Antibody Item No. 120560

Overview and Properties

Contents:	This vial contains peptide affinity-purified polyclonal antibody.
Synonym:	Cysteinyl-Leukotriene Receptor 2 (N-Term)
Immunogen:	Peptide corresponding to the N-terminal region of human CysLT ₂
Cross Reactivity:	(+) CysLT ₂ ; (-) 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), and Western blot (WB); the recommended starting dilution for ELISA is 1:200, ~1:1,000 for FC, and 1:250 for WB. 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, 08/04/2021

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

Cysteinyl leukotriene receptor 2 (CysLT₂) is a G protein-coupled receptor that mediates the actions of leukotriene C₄ (LTC₄), LTD₄, and, to a lesser extent, LTE₄.¹ It is expressed primarily in the heart, placenta, spleen, lymph nodes, and peripheral blood leukocytes.^{1,2} Signaling induced by LTC₄ through the CysLT₂ receptor induces the production of chemokines via NF-κB and AP-1 activation in endothelial cells, indicating a role in inflammation.³ A CysLT₂ receptor containing a leucine-to-glutamine substitution at position 129 (L129Q) has been found in human uveal melanoma cells and is associated with tumorigenesis *in vivo*.^{4,5} SNPs in *CYSLTR2* are associated with aspirin intolerance in patients with asthma.⁶ Cayman's CysLT₂ Receptor (N-Term) Polyclonal Antibody can be used for ELISA, flow cytometry (FC), and Western blot (WB) applications.

References

1. 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).
2. Takasaki, J., Kamohara, M., Matsumoto, M., *et al.* The molecular characterization and tissue distribution of the human cysteinyl leukotriene CysLT₂ receptor. *Biochem. Biophys. Res. Commun.* **274**, 316-322 (2000).
3. Thompson, C., Cloutier, A., Bossé, Y., *et al.* Signaling by the cysteinyl-leukotriene receptor 2. Involvement in chemokine gene transcription. *J. Biol. Chem.* **283**(4), 1974-1984 (2008).
4. Moore, A.R., Ceraudo, E., Sher, J.J., *et al.* Recurrent activating mutations of G-protein-coupled receptor *CYSLTR2* in uveal melanoma. *Nat. Genet.* **48**(6), 675-680 (2016).
5. Ceraudo, E., Horioka, M., Mattheisen, J.M., *et al.* Direct evidence that the GPCR *CysLTR2* mutant causative of uveal melanoma is constitutively active with highly biased signaling. *J. Biol. Chem.* **296**, 100163 (2021).
6. Park, J.S., Chang, H.S., Park, C.-S., *et al.* Association analysis of cysteinyl-leukotriene receptor 2 (*CYSLTR2*) polymorphisms with aspirin intolerance in asthmatics. *Pharmacogenet. Genomics* **15**(7), 483-492 (2005).

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