

# PRODUCT INFORMATION

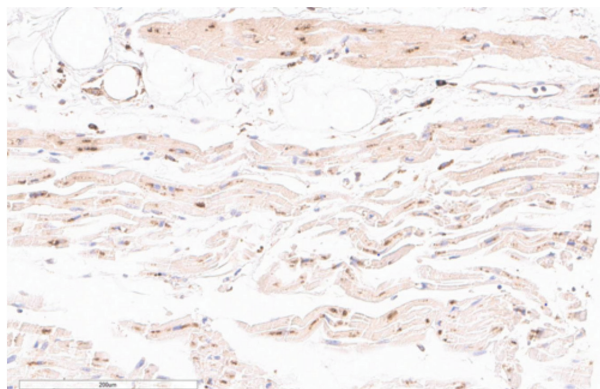


## CysLT<sub>2</sub> Receptor (N-Term) Polyclonal Antibody Item No. 120560

### Overview and Properties

<b>Contents:</b>	This vial contains peptide affinity-purified polyclonal antibody.
<b>Synonym:</b>	Cysteinyl-Leukotriene Receptor 2 (N-Term)
<b>Immunogen:</b>	Peptide corresponding to the N-terminal region of human CysLT <sub>2</sub>
<b>Cross Reactivity:</b>	(+) CysLT <sub>2</sub> ; (-) CysLT <sub>1</sub>
<b>Species Reactivity:</b>	(+) Human; (-) Rat
<b>Uniprot No.:</b>	Q9NS75
<b>Form:</b>	Lyophilized
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	TBS, pH 7.4, when reconstituted in 500 µl deionized water
<b>Host:</b>	Rabbit
<b>Applications:</b>	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 CysLT<sub>2</sub> 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, 11/07/2023

**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

# PRODUCT INFORMATION



## Description

---

Cysteinyl leukotriene receptor 2 (CysLT<sub>2</sub>) is a G protein-coupled receptor that mediates the actions of leukotriene C<sub>4</sub> (LTC<sub>4</sub>), LTD<sub>4</sub>, and, to a lesser extent, LTE<sub>4</sub>.<sup>1</sup> It is expressed primarily in the heart, placenta, spleen, lymph nodes, and peripheral blood leukocytes.<sup>1,2</sup> Signaling induced by LTC<sub>4</sub> through the CysLT<sub>2</sub> receptor induces the production of chemokines via NF-κB and AP-1 activation in endothelial cells, indicating a role in inflammation.<sup>3</sup> A CysLT<sub>2</sub> 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*.<sup>4,5</sup> SNPs in *CYSLTR2* are associated with aspirin intolerance in patients with asthma.<sup>6</sup> Cayman's CysLT<sub>2</sub> 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 CysLT<sub>2</sub> 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