

# PRODUCT INFORMATION

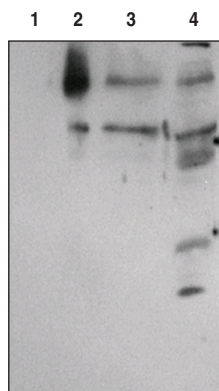


## BLT<sub>1</sub> Receptor Polyclonal Antibody Item No. 120114

### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
<b>Synonyms:</b>	BLTR <sub>1</sub> , Leukotriene B <sub>4</sub> Receptor 1, LTB <sub>4</sub> Receptor 1
<b>Immunogen:</b>	Synthetic peptide from the C-terminal region of human BLT <sub>1</sub>
<b>Cross Reactivity:</b>	(-) BLT <sub>2</sub> , CysLT <sub>1</sub> , and CysLT <sub>2</sub> Receptors 1
<b>Species Reactivity:</b>	(+) Human, mouse, and bovine BLT <sub>1</sub>
<b>Uniprot No.:</b>	Q15722
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	Flow cytometry (FC), Immunocytochemistry (ICC), Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



**Lane 1:** CHO lysate (30 µg)  
**Lane 2:** LTB<sub>4</sub> receptor transfected CHO cell lysate (30 µg)  
**Lane 3:** U937 cell lysate (50 µg)  
**Lane 4:** Bovine lung membrane (100 µg)

**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

---

The leukotriene B<sub>4</sub> receptor 1 (BLT<sub>1</sub> receptor), cloned from HL-60 human leukemia cells, has 352 amino acids and seven putative membrane-spanning domains.<sup>1</sup> The primary structure of the receptor is identical to that of a putative purinoceptor, P2Y<sub>7</sub>, which binds to micromolar concentrations of ATP.<sup>2</sup> Northern blotting reveals that the BLT<sub>1</sub> receptor is highly expressed in leukocytes, U937 cells, and to a much lower extent in spleen and thymus.<sup>1</sup> Sheep lung membranes have also been identified as a rich source for receptor isolation and purification.<sup>3</sup> A second LTB<sub>4</sub> receptor, BLT<sub>2</sub>, has recently been cloned and characterized.<sup>4-6</sup> This antibody was made against a peptide from the C-terminus of the BLT<sub>1</sub> receptor, which is located on the intracellular side of the plasma membrane. Therefore, when performing studies on whole cells, permeabilization of the cells is required for the antibody to enter the cytosol.

## References

---

1. Yokomizo, T., Izumi, T., Chang, K., *et al.* A G-protein-coupled receptor for leukotriene B<sub>4</sub> that mediates chemotaxis. *Nature* **387(6633)**, 620-624 (1997).
2. Akbar, G.K.M., Dasari, V.R., Webb, T.E., *et al.* Molecular cloning of a novel P2 purinoceptor from human erythroleukemia cells. *J. Biol. Chem.* **271(31)**, 18363-18367 (1996).
3. Votta, B., Keefer, J., and Mong, S. Characterization of the soluble leukotriene B<sub>4</sub> receptor from sheep lung membranes. *Biochem. J.* **270(1)**, 213-218 (1990).
4. Yokomizo, T., Kato, K., Terawaki, K., *et al.* A second leukotriene B<sub>4</sub> receptor, BLT2: A new therapeutic target in inflammation and immunological disorders. *J. Exp. Med.* **193(3)**, 421-431 (2000).
5. Kamohara, M., Takasaki, J., Matsumoto, M., *et al.* Molecular cloning and characterization of another leukotriene B<sub>4</sub> receptor. *J. Biol. Chem.* **275(35)**, 27000-27004 (2000).
6. Wang, S., Gustafson, E., Pang, L., *et al.* A novel hepatointestinal leukotriene B<sub>4</sub> receptor. *J. Biol. Chem.* **275(52)**, 40686-40694 (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