

# PRODUCT INFORMATION

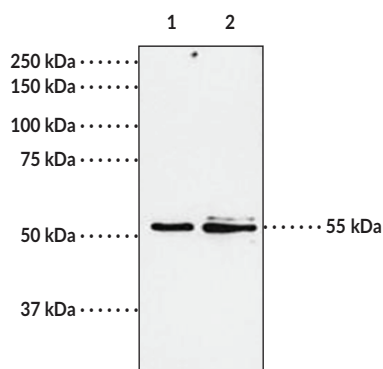


## BLT<sub>1</sub> Receptor Polyclonal Antiserum Item No. 100019

### Overview and Properties

<b>Contents:</b>	This vial contains polyclonal antiserum.
<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>Species Reactivity:</b>	(+) Human and bovine; (-) Mouse
<b>Uniprot No.:</b>	Q15722
<b>Form:</b>	Lyophilized
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥2 years
<b>Storage Buffer:</b>	Polyclonal antiserum when reconstituted in 100 µl of deionized water
<b>Host:</b>	Rabbit
<b>Applications:</b>	Confocal microscopy analysis (CF), flow cytometry (FC), immunohistochemistry (IHC) (formalin-fixed paraffin-embedded tissue), and Western blot (WB); the recommended starting dilution for CF and FC is 1:2,000, 1:200 for IHC (formalin-fixed paraffin-embedded tissue), and 1:1,000 for WB (80 and 60 kDa band pattern). <sup>2</sup> The BLT <sub>1</sub> receptor blocking peptide (Item No. 120112) is available for negative control experiments.

### Image



Lane 1: Human-kidney supernatant (25 µg)  
Lane 2: Human-kidney supernatant (50 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, 10/03/2018

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>3</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>4</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>3</sup> Sheep lung membranes have also been identified as a rich source for receptor isolation and purification.<sup>5</sup> A second LTB<sub>4</sub> receptor, BLT<sub>2</sub>, has recently been cloned and characterized.<sup>6-8</sup>

Cayman's BLT<sub>1</sub> receptor polyclonal antiserum is 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. Boie, Y., Stocco, R., Sawyer, N., *et al.* *Eur. J. Pharmacol.* **380**, 203-213 (1999).
2. Hennig, R., Ding, X.-Z., Tong, W.-G., *et al.* *Amer. J. Pathol.* **161**(2), 421-428 (2002).
3. Yokomizo, T., Izumi, T., Chang, K., *et al.* *Nature* **387**, 620-624 (1997).
4. Akbar, G.K.M., Dasari, V.R., Webb, T.E., *et al.* *J. Biol. Chem.* **271**, 18363-18367 (1996).
5. Votta, B., Keefer, J., and Mong, S. *Biochem. J.* **270**, 213-218 (1990).
6. Yokomizo, T., Kato, K., Terawaki, K., *et al.* *J. Exp. Med.* **193**, 421-431 (2000).
7. Kamohara, M., Takasaki, J., Matsumoto, M., *et al.* *J. Biol. Chem.* **275**, 27000-27004 (2000).
8. Wang, S., Gustafson, E., Pang, L., *et al.* *J. Biol. Chem.* **275**, 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