

# Product Information



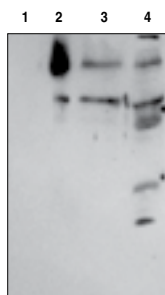
## BLT<sub>1</sub> Receptor Polyclonal Antibody

Item No. 120114 • Lot. No. XXXXX

<b>Synonyms:</b>	BLTR <sub>1</sub> ; Leukotriene B <sub>4</sub> Receptor 1; LTB <sub>4</sub> Receptor 1
<b>Contents:</b>	This vial contains (100-500 µg of peptide affinity-purified IgG, <i>lot specific</i> ) in 500 µl TBS, pH <i>lot specific</i> , containing <i>lot specific</i> % BSA and 0.02% sodium azide.
<b>Host:</b>	Rabbit
<b>Antigen:</b>	Human BLT <sub>1</sub> receptor amino acids 331-352 (ALEPGPSESILTASSPLKLNELN) <sup>1</sup>
<b>Cross-reactivity:</b>	(+) Human and bovine BLT <sub>1</sub> receptor; (-) Mouse BLT <sub>1</sub> receptor. Homology of peptide antigen with BLT <sub>1</sub> receptor from other species: Human: A L E P G P S E S I L T A S S P L K L N E L N Mouse: c p E P G P t d S f m t S S t i p e s - s k Guinea pig: A p E P G a S g S L d g - - L K q s E s d
<b>Stability:</b>	≥ 1 year at -20°C
<b>Applications:</b>	Use at a dilution of 1: <i>lot specific</i> for western blot or 1: <i>lot specific</i> for flow cytometry. The antibody can also be used for immunocytochemistry and immunohistochemistry.
<b>Concentration:</b>	Varies by lot, from 0.2-1.0 mg/ml (100-500 µg/vial). Always 100 µl final working volume for western blotting.

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. For flow cytometry and confocal microscopy analysis of the receptor, the antibody can be diluted approximately 1: *lot specific*; use at a dilution of approximately 1: *lot specific* for western blot analysis.



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)

### References

1. Yokomizo, T., Izumi, T., Chang, K., *et al. Nature* **387**, 620-624 (1997).
2. Akbar, G.K.M., Dasari, V.R., Webb, T.E., *et al. J. Biol. Chem.* **271**, 18363-18367 (1996).
3. Votta, B., Keefer, J., Mong, S. *Biochem. J.* **270**, 213-218 (1990).
4. Yokomizo, T., Kato, K., Terawaki, K., *et al. J. Exp. Med.* **193**, 421-431 (2000).
5. Kamohara, M., Takasaki, J., Matsumoto, M., *et al. J. Biol. Chem.* **275**, 27000-27004 (2000).
6. Wang, S., Gustafson, E., Pang, L., *et al. J. Biol. Chem.* **275**, 40686-40694 (2000).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/120114](http://www.caymanchem.com/catalog/120114)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/20/2011

## Cayman Chemical

### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

### Phone

(800) 364-9897  
(734) 971-3335

### Fax

(734) 971-3640

### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

### Web

[www.caymanchem.com](http://www.caymanchem.com)