

# Product Information



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

Catalog No. 120111 • Lot. No. XXXX

<b>Contents:</b>	This vial contains XXX µg of protein-A purified IgG in XXX µl of PBS, pH 7.4, containing 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide
<b>Synonyms:</b>	BLTR <sub>1</sub> ; Leukotriene B <sub>4</sub> Receptor 1; LTB <sub>4</sub> Receptor 1
<b>Antigen:</b>	HeLa cells transfected with human BLT <sub>1</sub>
<b>Host:</b>	mouse, clone 7B1
<b>Cross Reactivity:</b>	(+) Human BLT <sub>1</sub> receptor; (-) BLT <sub>2</sub> , CysLT <sub>1</sub> , and CysLT <sub>2</sub> receptors <sup>1</sup>
<b>Stability:</b>	≥1 year at -20°C
<b>Applications:</b>	Flow cytometry, IHC of frozen sections, and ICC at XX µg/ml. BLT <sub>1</sub> antagonism at XX µg/ml with semi-purified receptor preparations. The best working dilution of the antibody may need to be determined empirically for each application. Does not work for western blotting.

The BLT<sub>1</sub> receptor, cloned from HL-60 human leukemia cells, has 352 amino acids and seven putative membrane-spanning domains.<sup>2</sup> The primary structure of the receptor is identical to that of a putative purinoceptor, P2Y7, which binds to micromolar concentrations of ATP.<sup>3</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>2</sup> Sheep lung membranes have also been identified as a rich source for receptor isolation and purification.<sup>4</sup> A second LTB<sub>4</sub> receptor, BLT<sub>2</sub>, has also been cloned and characterized.<sup>5-7</sup>

### References

- Pettersson, A., Boketoft, A., Sabirsh, A., *et al.* First-generation monoclonal antibodies identifying the human leukotriene B<sub>4</sub> receptor-1. *Biochem. Biophys. Res. Commun.* **279**, 520-525 (2000).
- Yokomizo, T., Izumi, T., Chang, K., *et al.* A G-protein-coupled receptor for leukotriene B<sub>4</sub> that mediates chemotaxis. *Nature* **387**, 620-624 (1997).
- Akbar, G.K.M., Dasari, V.R., Webb, T.E., *et al.* Molecular cloning of a novel P<sub>2</sub> purinoceptor from human erythroleukemia cells. *J. Biol. Chem.* **271**, 18363-18367 (1996).
- Votta, B., Keefer, J., and Mong, S. Characterization of the soluble leukotriene B<sub>4</sub> receptor from sheep lung membranes. *Biochem. J.* **270**, 213-218 (1990).
- Yokomizo, T., Kato, K., Terawaki, K., *et al.* A second leukotriene B<sub>4</sub> receptor, BLT<sub>2</sub>: A new therapeutic target in inflammation and immunological disorders. *J. Exp. Med.* **193**, 421-431 (2000).
- Kamohara, M., Takasaki, J., Matsumoto, M., *et al.* Molecular cloning and characterization of another leukotriene B<sub>4</sub> receptor. *J. Biol. Chem.* **275**, 27000-27004 (2000).
- Wang, S., Gustafson, E., Pang, L., *et al.* A novel hepatointestinal leukotriene B<sub>4</sub> receptor. *J. Biol. Chem.* **275**, 40686-40694 (2000).

### Related Products

BLT<sub>1</sub> Receptor Polyclonal Antiserum - Cat. No. 100019 • BLT<sub>1</sub> Receptor Polyclonal Antibody - Cat. No. 120114 • BLT<sub>2</sub> Receptor Polyclonal Antibody - Cat. No. 120124

### 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

**Web**  
www.caymanchem.com

**WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.**

#### 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 under separate cover to the MSDS supervisor at 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, 06/27/2006