

# PRODUCT INFORMATION



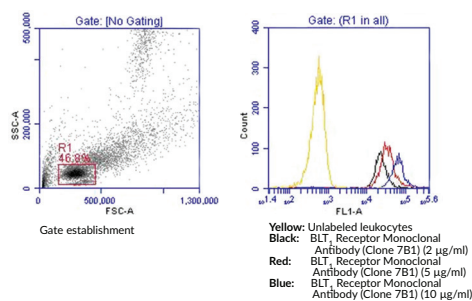
## BLT<sub>1</sub> Receptor Monoclonal Antibody (Clone 7B1)

Item No. 120111

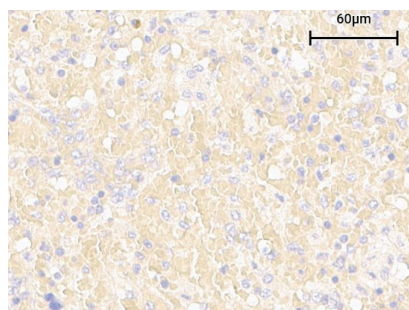
### Overview and Properties

<b>Contents:</b>	This vial contains 200 µg of protein G-purified monoclonal antibody.
<b>Synonyms:</b>	BLTR <sub>1</sub> , Leukotriene B <sub>4</sub> Receptor 1, LTB <sub>4</sub> Receptor 1
<b>Immunogen:</b>	HeLa cells transfected with human BLT <sub>1</sub>
<b>Cross Reactivity:</b>	(+) BLT <sub>1</sub> , (-) BLT <sub>2</sub> , CysLT <sub>1</sub> , and CysLT <sub>2</sub> receptors 1 <sup>1</sup>
<b>Species Reactivity:</b>	(+) Human
<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, 0.1% BSA, and 0.02% sodium azide
<b>Clone:</b>	7B1
<b>Host:</b>	Mouse
<b>Isotype:</b>	IgG2a
<b>Applications:</b>	Flow cytometry (FC), Immunofluorescence (IF), and Immunohistochemistry (IHC). BLT <sub>1</sub> antagonism at 5 µg/ml with semi-purified receptor preparations; the recommended starting concentration is 5 µg/ml for FC and IHC; the recommended starting dilution is 5 µg/ml for IF. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. Does not work for Western blotting.

### Images



Whole human blood was fixed with 4% formaldehyde. Erythrocytes were lysed by the addition of 0.1% Triton-X. Cells were blocked with 3% normal goat serum then treated sequentially with primary (BLT<sub>1</sub> Receptor Monoclonal Antibody (Clone 7B1)) and secondary (goat anti-mouse FITC labeled) antibodies with 1 hour, room temperature incubations. The samples were analyzed using an Accuri flow cytometer using the FL1 detector. Color compensation: FL2 by FL1: 5.5.



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human spleen tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with BLT<sub>1</sub> Monoclonal Antibody (Clone 7B1) (Item No. 120111) at a 1:80 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, 11/08/2023

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 human BLT<sub>1</sub> receptor is a G protein-coupled receptor that mediates the proinflammatory effects of leukotriene B<sub>4</sub> (LTB<sub>4</sub>).<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>2</sup> Sheep lung membranes have also been identified as a rich source for receptor isolation and purification.<sup>3</sup> A second, low-affinity LTB<sub>4</sub> receptor, BLT<sub>2</sub>, has also been cloned and characterized.<sup>4-6</sup> Cayman Chemical's BLT<sub>1</sub> Receptor Monoclonal Antibody is a useful tool for the detection of human BLT1 by flow cytometry, immunofluorescence, and immunohistochemistry. The antibody does not cross react with the other leukotriene receptors (BLT<sub>2</sub>, CysLT<sub>1</sub>, or CysLT<sub>2</sub>) and does not work for Western blot analysis of BLT<sub>1</sub>.

## References

---

1. 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(2)**, 520-525 (2000).
2. 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).
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