

Product Information



BLT₂ Receptor Polyclonal Antibody

Item No. 120124 • Lot. No. XXXXX

Synonyms:	BLTR ₂ ; Leukotriene B ₄ Receptor 2; LTB ₄ Receptor 2
Contents:	This vial contains (100-500 µg of peptide affinity-purified IgG, <i>lot specific</i>) in 500 µl TBS, pH 7.4, containing <i>lot specific</i> mg/ml BSA, <i>lot specific</i> % glycerol, and <i>lot specific</i> % sodium azide.
Host:	Rabbit
Antigen:	The BLT ₂ receptor polyclonal antibody was raised against a synthetic peptide from the human BLT ₂ receptor amino acids 325-338 (MELRTTPQLKVVGQ) conjugated to KLH. ¹⁻³
Cross Reactivity:	(+) Human and mouse BLT ₂ receptor; other species not tested. Human MELRTTPQLKVVGQ Mouse MELRTTP k LKVMGQ
Stability:	≥2 years at -20°C
Application:	Recommended starting dilutions for western blotting: <i>lot specific:lot specific</i> (<i>lot specific</i> µg/ml) (recommended primarily for transfected cells). ⁴
Concentration:	Varies by lot, from 0.2-1.0 mg/ml (100-500 µg/vial). Always 100 ml final working volume for western blotting.

Leukotriene B₄ (LTB₄) promotes a number of leukocyte functions including aggregation, stimulation of ion fluxes, superoxide anion production, chemotaxis, and chemokinesis. Two G-protein coupled receptors, BLT₁ and BLT₂, which mediate the actions of LTB₄ have been cloned and characterized.^{1-3,5} The BLT₁ receptor is expressed in leukocytes, U937 cells, and to a much lower extent in spleen and thymus.⁵ The human BLT₂ receptor is a 358 amino acid protein with approximately 40% homology to BLT₁ at the amino acid level.¹⁻³ The BLT₂ receptor is more broadly expressed than the BLT₁ receptor, with highest levels observed in liver, spleen, ovary, intestine, and peripheral blood leukocytes.¹⁻³ Activation of the BLT₂ receptor by LTB₄ leads to Ca²⁺ mobilization as well as decreases in cyclic AMP, indicating interaction of the receptor with G_q and G_i classes of G-proteins.²



References

1. Yokomizo, T., Kato, K., Terawaki, K., *et al.* A second leukotriene B₄ receptor, BLT₂: A new therapeutic target in inflammation and immunological disorders. *J. Exp. Med.* **193**, 421-431 (2000).
2. Wang, S., Gustafson, E., Pang, L., *et al.* A novel hepatointestinal leukotriene B₄ receptor. *J. Biol. Chem.* **275**, 40686-40694 (2000).
3. Kamohara, M., Takasaki, J., Matsumoto, M., *et al.* Molecular cloning and characterization of another leukotriene B₄ receptor. *J. Biol. Chem.* **275**, 27000-27004 (2000).
4. Paredes, Y., Massicotte, F., Pelletier, J.-P., *et al.* Study of the role of leukotriene B₄ in abnormal function of human subchondral osteoarthritis osteoblasts. Effects of cyclooxygenase and/or 5-lipoxygenase inhibition. *Arthritis Rheuma (Munch)* **46(7)**, 1804-1812 (2002).
5. Yokomizo, T., Izumi, T., Chang, K., *et al.* A G-protein-coupled receptor for leukotriene B₄ that mediates chemotaxis. *Nature* **387**, 620-624 (1997).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/120124

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

Web

www.caymanchem.com