

Product Information



Toll-Like Receptor 4 Monoclonal Antibody (Clone HTA125)

Catalog No. 13589

Contents:	This vial contains 100 µg of protein G-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
Synonym:	TLR4
Antigen:	Ba/F3 cell line expressing human TLR4 cell surface antigen
Isotype:	IgG _{2ak}
Host:	Mouse, clone HTA125
Cross Reactivity:	(+) Canine and human TLR4
Storage:	≤6 months at 4°C; ≥6 months at -20°C
Applications:	Flow cytometry (FC) (intracellular and cell surface), immunoprecipitation (IP), neutralization, and immunocytochemistry. The recommended starting concentration for FC (cell surface) is 2-3 µg/1x10 ⁶ cells.

The toll-like receptors (TLRs) in mammals comprise a family of transmembrane proteins characterized by multiple copies of leucine rich repeats in the extracellular domain and an interleukin-1 (IL-1) receptor motif in the cytoplasmic domain. Like their counterparts in *Drosophila*, TLRs signal through adaptor molecules.¹ The TLR family is a phylogenetically conserved mediator of innate immunity that is essential for microbial recognition.² Most mammalian species have between ten and fifteen types of TLRs. Ten functional TLRs (TLR1-10) have been identified in human. Humans also encode a TLR11 gene but it contains several stop codons and protein is not expressed. However, mouse and rat TLR11 are functional, and it is thought that human TLR11 function was lost during evolution. Historically speaking, TLR expression has been most extensively studied in the immune system. Overall, TLRs are highly expressed in immune competent cells, including macrophages, dendritic cells, neutrophils, mucosal epithelial cells and dermal endothelial cells. However, TLRs have also been identified in many other cell types and anatomical tissue locations where they are expressed either constitutively or induced during infection. Among this family of receptors TLR2 and TLR4 have been most studied. These studies have suggested that TLR2 and TLR4 may serve as potential main mediators of LPS signaling.^{3,4} The TLR4 cDNA codes for a protein consisting of 799 amino acids with approximate molecular weight of 88 kDa.⁵

References

1. Srivastava, M.D., Srivastava, B.I.S. Expression of mRNA and proteins for toll-like receptors, associated molecules, defensins and LL-37 by SRIK-NKL, and CD⁸ NK/T cell line. *Leukemia Research* **29**(7), 813-820 (2005).
2. Gibson, F.C., Hong, C., Chou, H.-H., *et al.* Innate immune recognition of invasive bacteria accelerates atherosclerosis in apolipoprotein E-deficient mice. *Circulation* **109**, 2801-2806 (2004).
3. Cognasse, F., Hamzeh, H., Chavarin, P., *et al.* Evidence of toll-like receptor molecules on human platelets. *Immunology & Cell Biology* **83**, 196-198 (2005).
4. Wong, C.K., Cheung, P.F.Y., Ip, W.K., *et al.* Intracellular signaling mechanisms regulating toll-like receptor-mediated activation of eosinophils. *Am. J. Respir. Cell and Mol. Biol.* **37**, 85-96 (2007).
5. Scheel, O., Papavlassopoulos, M., Clunck, R., *et al.* Cell activation by ligands of the toll-like receptor and interleukin-1-receptor family depends on the function of the large-conductance potassium channel maxiK in human macrophages. *Infection and Immunity* **74**(7), 4354-4356 (2006).

Related Products

Toll-Like Receptor 1 Polyclonal Antibody - Cat. No. 13582 • Toll-Like Receptor 2 Monoclonal Antibody (Clone TL2.1) - Cat. No. 13587 • Toll-Like Receptor 3 Monoclonal Antibody (Clone 40C1285.6) - Cat. No. 13588 • Toll-Like Receptor 5 Monoclonal Antibody (Clone 85B152.5) - Cat. No. 13595 • Toll-Like Receptor 6 Polyclonal Antibody - Cat. No. 13590 • Toll-Like Receptor 7 Polyclonal Antibody - Cat. No. 13591 • Toll-Like Receptor 8 Monoclonal Antibody (Clone 44C143) - Cat. No. 13592 • Toll-Like Receptor 9 Monoclonal Antibody (Clone 26C593.2) - Cat. No. 13593 • Toll-Like Receptor 10 Monoclonal Antibody - Cat. No. 13584 • Toll-Like Receptor 11 Polyclonal Antibody - Cat. No. 13585 • Toll-Like Receptor 12 Polyclonal Antibody - Cat. No. 13586

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, 12/01/2009

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