

Product Information



MyD88 Polyclonal Antibody

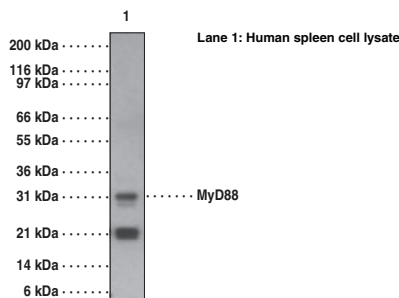
Item No. 13746

- Contents:** This vial contains 100 µg of protein G-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
- Antigen:** Synthetic peptide corresponding to human MyD88 amino acids 233-248
- Host:** Rabbit
- Cross Reactivity:** (+) Human, mouse, and rat MyD88
- Storage:** ≤6 months at 4°C; ≥6 months at -20°C
- Application:** Western blot (WB). The recommended starting concentration for WB is 0.5-2 µg/ml.

MyD88, a protein involved in interleukin-1 (IL-1) mediated signaling, was originally isolated as myeloid differentiation primary response gene.^{1,2} MyD88 possesses an N-terminal death domain similar to cytoplasmic segments of TNF receptor 1, Fas, and a C-terminal region related to IL-1 and toll receptors. Overexpression of MyD88 induces activation of the c-Jun N-terminal kinase and NF-κB through its death domain.^{3,4} The C-terminus of MyD88 interacts with the IL-1 receptor and blocks NF-κB activation induced by IL-1, but not by TNF.

References

1. Bonnert, T.P., Garka, K.E., Parnet, P., *et al.* The cloning and characterization of human MyD88: A member of an IL-1 receptor related family. *FEBS Lett.* **402(1)**, 81-84 (1997).
2. Hardiman, G., Rock, F.L., Balsubramanian, S., *et al.* Molecular characterization and modular analysis of human MyD88. *Oncogene* **13(11)**, 2467-2475 (1996).
3. Burns, K., Martinon, F., Esslinger, C., Pahl, H., *et al.* MyD88, an adapter protein involved in interleukin-1 signaling. *J.Biol.Chem.* **273(20)**, 12203-12209 (1998).
4. Wesche, H., Henzel, W.J., Shillinglaw, W., Li, S., *et al.* MyD88: An adapter that recruits IRAK to the IL-1 receptor complex. *Immunity* **7(6)**, 837-847 (1997).



Related Products

Interleukin-1α (human) EIA Kit - Item No. 583301 • Interleukin-1β (human) EIA Kit - Item No. 583311

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 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, 07/07/2010