

Product Information

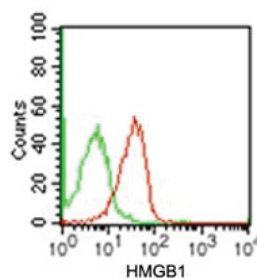


HMGB1 Monoclonal Antibody (Clone IMG19N15F4)

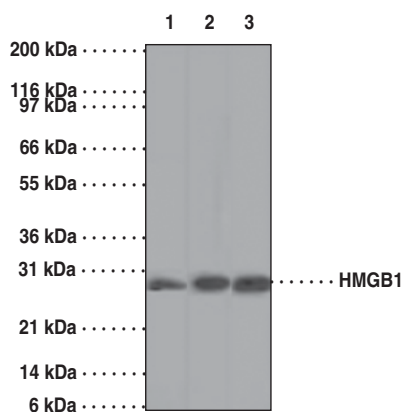
Item No. 11513

- Contents:** This vial contains 100 µg of Protein G purified IgG_{1κ} in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
- Synonyms:** High Mobility Group Protein B1, HMG1
- Antigen:** Full length recombinant human HMGB1
- Host:** Mouse, clone IMG19N15F4
- Isotype:** IgG_{1κ}
- Cross Reactivity:** (+) Human and mouse HMGB1; other species not tested.
- Stability:** ≥1 year at -20°C
- Applications:** Flow cytometry (FC), immunohistochemistry (IHC) (paraffin-embedded tissue), and western blot (WB); the recommended starting concentration for FC is 0.5 µg/5x10⁵ cells, 5 µg/ml for IHC, and 1-3 µg/ml for WB. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

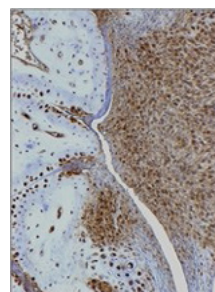
High Mobility Group Protein B1 (HMGB1) belongs to the HMGB family and contains two HMG box DNA-binding domains. It is a highly conserved, ubiquitous protein present in the nuclei and cytoplasm of nearly all cell types, and is a necessary and sufficient mediator of inflammation during sterile and infection-associated responses. HMGB1 also acts as a DNA nuclear binding protein that has been shown to be an early trigger of sterile inflammation in animal models of trauma-hemorrhage *via* the activation of the Toll-like receptor 4 (TLR4) and the receptor for the advanced glycation endproducts (RAGE).¹ It is reported that the level of HMGB1 is elevated during sterile tissue injury, infection, lethal endotoxemia or sepsis, collagen-induced arthritis, and ischemia-reperfusion induced tissue injury.²



Intracellular flow cytometric analysis using HMGB1 antibody. Human Jurkat cells were probed using 0.5 µg of HMGB1 antibody (red) and 0.5 µg of isotype control antibody (green).



Western blot analysis using HMGB1 antibody at 2 µg/ml.
Lane 1: Full-length human HMGB1
Lane 2: Human Jurkat cell lysate
Lane 3: Mouse NIH 3T3 cell lysate



Immunohistochemical analysis of formalin-fixed, decalcified, paraffin embedded tissue sections from the paws of mice with collagen-induced arthritis using the HMGB1 antibody. The antibody was used at 5 µg/ml. Data courtesy of Dr. Ulf Andersson, Karolinska Institute.

References

1. Cohen, M.J., Brohi, K., Calfee, C.S., *et al.* *Crit. Care* **13**(6), 1-10 (2009).
2. Yang, H. and Tracey, K.J. *Biochim. Biophys. Acta* **1799**, 149-156 (2010).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/11513

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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 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, 01/14/2014

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