

PRODUCT INFORMATION



HMGB1 Polyclonal Antibody (aa 100-150)

Item No. 11515

Overview

Contents: This vial contains 100 µg of Protein A purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.

Synonyms: High Mobility Group Protein B1, HMG1

Immunogen: Human HMGB1 amino acids 100-150

Species Reactivity: (+) Human, bovine, chicken, mouse, New World monkey, and rat HMGB1; other species not tested.

Properties

Storage: -20°C (as supplied)

Stability: ≥1 year

Host: Rabbit

Applications

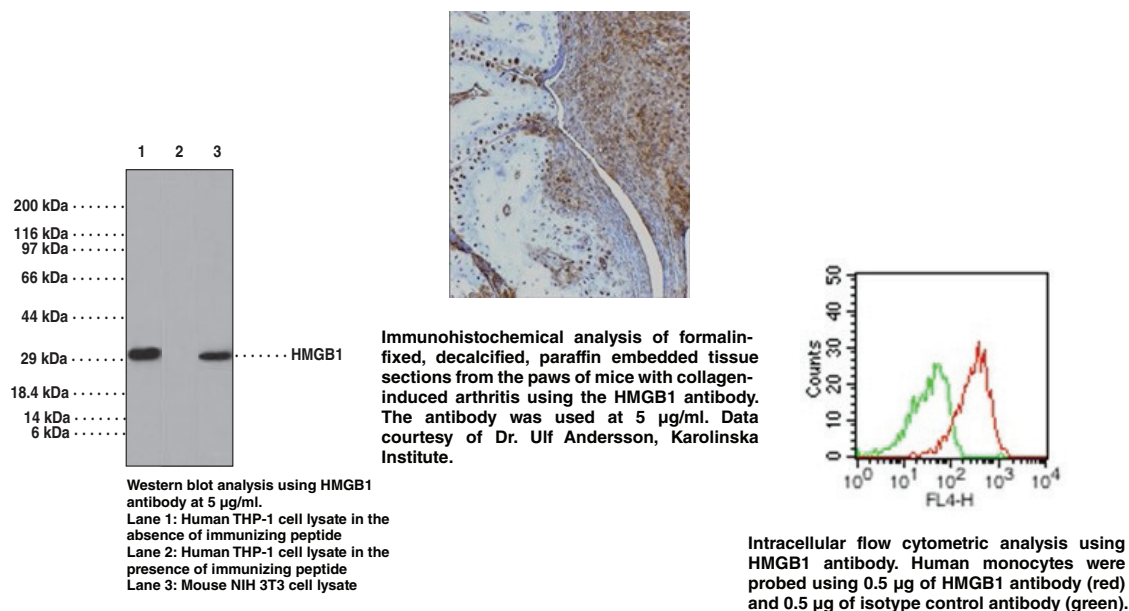
Flow cytometry: Suggested starting concentration is 0.5 µg/10⁶ cells

Immunohistochemistry (paraffin-embedded tissue): Suggested starting concentration is 5 µg/ml

Western blot: Suggested starting concentration is 5-10 µg/ml

Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Images



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, 07/30/2015

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