

HMGB1 (human) monoclonal antibody (Moby-2)

High mobility group 1 protein (HMGB1) is a chromatin component that is either passively released from necrotic cells or actively secreted in response to angiogenic and inflammatory signals. HMGB1 itself may signal through the receptor for advanced glycation end products (RAGE), and via the Toll-like receptors, TLR2 and TLR4. HMGB1 functions as a late mediator of immune response. Apoptotic cells also release HMGB1 but in an oxidized form that does not trigger the immune response.

This antibody is covered by our [Worry-Free Guarantee](#).

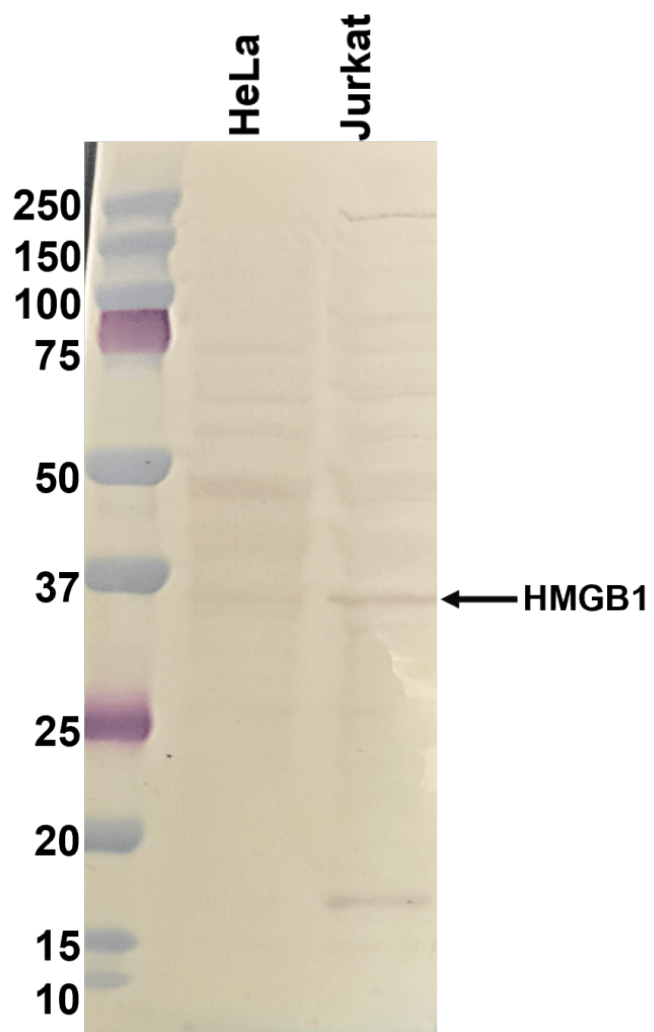
Ordering Information

[Order Online »](#)

ALX-804-873-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HMGB1 mAb (Prod. No. ALX-804-873) at 1:200 dilution. The HRP-Conjugated anti-mouse IgG was used as the secondary antibody. Detection was with colorimetric substrate.

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	High mobility group protein B1
Application	ELISA, IP, WB
Clone	Moby-2
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human HMGB1.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Immunoprecipitation (1:200)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Human
UniProt ID	P09429

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 8, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)