

# PRODUCT INFORMATION



## HMGB1 (human, recombinant)

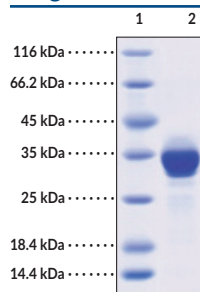
Item No. 43701

### Overview and Properties

<b>Synonyms:</b>	Amphoterin, High-mobility Group Box 1 Protein, High Mobility Group Protein B1, HMG-1, SBP-1
<b>Source:</b>	Active recombinant human C-terminal His-tagged HMGB1 expressed in HEK293 cells
<b>Amino Acids:</b>	2-215 (full length)
<b>Uniprot No.:</b>	P09429
<b>Molecular Weight:</b>	26.21 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Purity:</b>	≥95% estimated by SDS-PAGE
<b>Supplied in:</b>	Lyophilized from sterile PBS, pH 7.4
<b>Endotoxin Testing:</b>	<1.0 EU/μg, determined by the LAL endotoxin assay
<b>Bioactivity:</b>	See figures for details
<b>Specific Activity:</b>	Batch specific

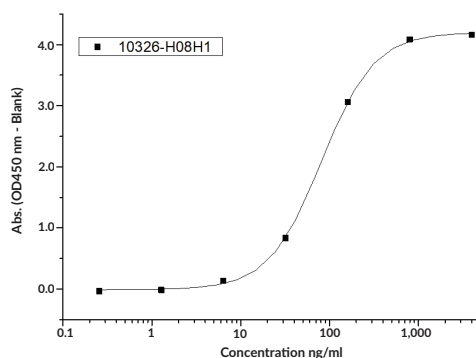
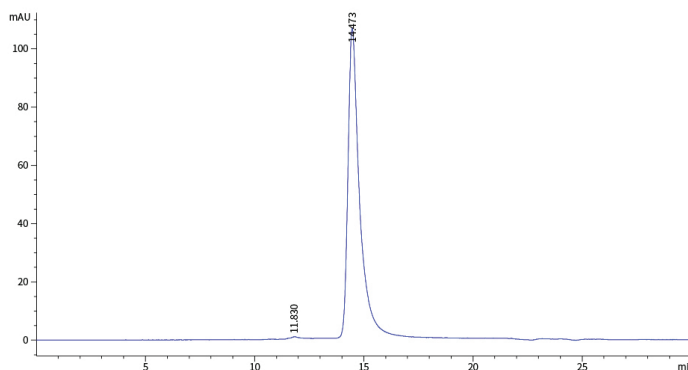
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Images



Lane 1: MW Markers  
Lane 2: HMGB1

**SDS-PAGE Analysis of HMGB1.** This protein has a calculated molecular weight of 26.21 kDa. It has an apparent molecular weight of approximately 33.2 kDa by SDS-PAGE under reducing conditions due to glycosylation.



Immobilized HMGB1 (human, recombinant) at 2 μg/ml (100 μl/well) can bind Recombinant Human AGER / RAGE Protein (Fc Tag), with an EC<sub>50</sub> value of 40-120 ng/ml.

**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, 12/12/2025

**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

# PRODUCT INFORMATION



## Description

---

High mobility group protein B1 (HMGB1) is a member of the HMGB subfamily of proteins with roles in promoting proper DNA structure and mediating extracellular inflammation.<sup>1,2</sup> It is composed of two HMG box DNA-binding domains and a negatively charged C-terminus.<sup>1,3</sup> HMGB1 is ubiquitously expressed and found in the nucleus and extracellular space.<sup>1</sup> In the nucleus, it binds DNA and histones and ensures proper DNA structure.<sup>3</sup> HMGB1 is secreted by cells in response to regulated cell death pathways, tissue damage, ischemia, organ transplants, or pathogen infection, where it further increases local inflammation and stimulates additional HMGB1 secretion by immune cells.<sup>1,2,4</sup> Serum levels of HMGB1 are increased in patients with systemic inflammatory response syndrome (SIRS), a condition characterized by immune system hyperreactivity to infectious and non-infectious stimuli, and a SNP in the promoter region of *HMGB1* is associated with poor survival in SIRS patients.<sup>5</sup> Cayman's HMGB1 (human, recombinant) protein can be used in binding assays. This protein consists of 225 amino acids and has a calculated molecular weight of 26.21 kDa. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is approximately 33.2 kDa due to glycosylation.

## References

---

1. Tang, D., Kang, R., Zeh, H.J., *et al.* The multifunctional protein HMGB1: 50 years of discovery. *Nat. Rev. Immunol.* **23**(12), 824-841 (2023).
2. Yang, H. and Tracey, K.J. Targeting HMGB1 in inflammation. *Biochim. Biophys. Acta* **1799**(1-2), 149-156 (2010).
3. Bianchi, M.E., Beltrame, M., and Paonessa, G. Specific recognition of cruciform DNA by nuclear protein HMG1. *Science* **243**(4894 Pt 1), 1056-1059 (1989).
4. Ku, S.K., Lee, I.C., Kim, J.A., *et al.* Anti-septic effects of pellicitorine in HMGB1-induced inflammatory responses *in vitro* and *in vivo*. *Inflammation* **37**(2), 338-348 (2014).
5. Kornblit, B., Munthe-Fog, L., Madsen, H.O., *et al.* Association of HMGB1 polymorphisms with outcome in patients with systemic inflammatory response syndrome. *Crit. Care* **12**(3), R83 (2008).

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