

# HMGB1 monoclonal antibody (KS1)

The chromatin non-histone high mobility group of proteins (HMG)-1 and -2 contain conserved domains known as HMG-boxes. Both HMG-1 and HMG-2 contain an N-terminal HMG box, a central HMG box and an acidic carboxy terminus. HMG-1/2 have been shown to facilitate the binding of various sequence-specific transcription factors to their respective DNA binding sites, such as steroid hormone receptors, and p53.

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

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	SBP-1, High-mobility group box 1
Application	WB
Application Notes	Detects a band of ~29kDa by Western blot.
Clone	KS1
Formulation	Liquid. In PBS containing 50% glycerol.
Host	Mouse
Immunogen	Pig HMGB1 (HMG1).
Isotype	IgG2a
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (5µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P12682
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>