

α -HMG1

anti-High Mobility Group 1
rabbit, polyclonal

Cat. No.	Amount
ABD-012	125 μ l

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied as high titer anti-serum.

Description

High mobility group 1 (HMG1) is a 26 kDa highly conserved non-sequence-specific DNA-binding nuclear protein. Mammalian HMG1 has two homologous DNA-binding domains HMG boxes, A and B (each of 80–90 amino-acid residues), linked by a short basic region to an acidic C-terminal domain containing 30 consecutive Asp and Glu residues. HMG1 has been implicated in a number of fundamental biological processes including transcription, replication and recombination, in which it plays a role in manipulating DNA structure by bending, looping, compaction or unwinding, or by direct contacts with distinct cellular proteins. HMG1 can act as a repressor, by interacting with TBP to block pre-initiation complex formation or as an activator, by facilitating the binding of various transcription factors to their cognate DNA sequences. Most recently, it was discovered that HMG1 is a late mediator of delayed endotoxin lethality by activating downstream cytokine release.

Specificity

α -HMG1 reacts with the HMG1 protein in HeLa nuclear extract (cat.# PR-777) by Western blotting. Recommended dilution range for Western blot analysis: 1:2000 - 1:5000.

Selected References:

- Bianchi et al. (1989) *Science* **243**:1056.
- Bustin et al. (1990) *Biochem. Biophys. Acta* **1049**:231.
- Zappavigna et al. (1996) *EMBO J.* **15**:4981.
- Ge et al. (1994) *J. Biol. Chem.* **269**:17136.
- Zlatanova et al. (1998) *FASEB J.* **12**:421.
- Stelzer et al. (1994) *Mol. Cell. Biol.* **14**:4712.
- Lu et al. (2000) *J. Biol. Chem.* **10**:35006.
- Jayaraman et al. (1998) *Genes Dev.* **12**:462.
- Onate et al. (1994) *Mol. Cell. Biol.* **14**:3376.
- Yang et al. (2001) *Shock* **15**:247.