

## Recombinant HMGB1 protein

---

**Catalog No:** 31247

**Expressed In:** *E. coli*

**Quantity:** 10 µg

**Concentration:** 0.15 µg/µl

**Source:** Human

**Buffer Contents:** 10 µg of recombinant protein supplied at a concentration of 0.15 µg/µl in 20 mM Tris-HCl, pH 7.3, 100 mM KCl, 0.2 mM EDTA and 20% glycerol.

**Background:** **HMGB1** (High Mobility Group Box 1) is a DNA binding proteins that associates with chromatin and has the ability to bend DNA. Binds preferentially single-stranded DNA. Involved in V(D)J recombination by acting as a cofactor of the RAG complex. Acts by stimulating cleavage and RAG protein binding at the 23 bp spacer of conserved recombination signal sequences (RSS). Heparin-binding protein that has a role in the extension of neurite-type cytoplasmic processes in developing cells (By similarity).

**Protein Details:** Recombinant HMGB1 protein was expressed in *E. coli* as the full-length human protein (accession number NP\_002119) with an N-terminal polyhistidine tag. The molecular weight of the protein is ~27 kDa.

**Application Notes:** Recombinant HMGB1 protein is suitable for use in *in vitro* transcription assays. We recommended to start with a range of protein concentrations (*e.g.* 5 ng, 50 ng, and 500 ng) to determine the optimal amount needed for each *in vitro* transcription assay.

**Storage and Guarantee:** Lyophilized proteins can be stored at -20°C or -80°C, preferably desiccated. Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.