

HMGN1 antibody (pAb)

Catalog Nos: 39833, 39834

RRID: AB_2793360

Isotype: IgG

Application(s): IP, WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 20 kDa

Background:

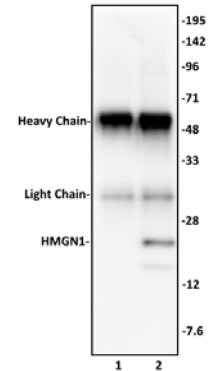
Immunogen: This HMGN1 antibody was raised against a peptide within the C-terminal region of human HMGN1.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

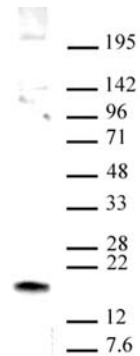
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

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



HMGN1 antibody (pAb) tested by Immunoprecipitation.

10 μ l of HMGN1 antibody was used to immunoprecipitate HMGN1 from 400 μ g of HeLa whole cell extract (lane 2). 10 μ l of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the HMGN1 antibody at a dilution of 1:500.



HMGN1 antibody (pAb) tested by Western blot.

Detection of HMGN1 by Western blot analysis. Whole cell extract of HeLa cells (20 μ g) probed with HMGN1 antibody at a dilution of 1:500.