

## HMGA2 / HMGI-C antibody (pAb)

**Catalog Nos:** 61041, 61042

**RRID:** AB\_2614960

**Isotype:** IgG

**Application(s):** IP, WB

**Reactivity:** Human, Mouse

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 20 kDa

**Background:** HMGA2 (High-mobility group AT-hook protein 2, HMGI-C) is a member of the HMG (High mobility group) protein family, which are non-histone chromatin-associated proteins involved in DNA organization and regulation of transcription. The HMG protein family consists of six proteins and is subdivided into three subfamilies: the HMG-1/-2 subfamily, the HMG-I/Y subfamily and the HMG-14/-17 subfamily. Mouse models of HMGA2 function indicate that it is involved in proliferation at the cellular level and obesity at the organismal level. Overexpression of HMGA2 is observed in a variety of cancers.

**Immunogen:** This HMGA2 antibody was raised against a peptide within the N-terminal region of human HMGA2 / HMGI-C.

**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:

Applications Validated by Active Motif:

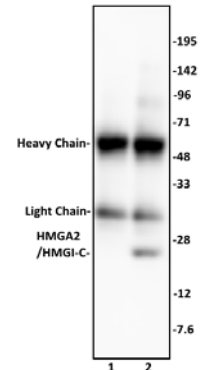
IP: 10 µl per IP

WB\*: 1:1,000 - 1:5,000 dilution

\*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

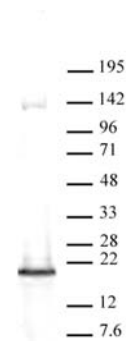
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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



### HMGA2 / HMGI-C antibody (pAb) tested by Immunoprecipitation.

10 µl of HMGA2 / HMGI-C antibody was used to immunoprecipitate HMGA2 / HMGI-C from 250 µg of HepG2 nuclear cell extract (lane 2). 10 µl of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the HMGA2 / HMGI-C antibody at a dilution of 1:3,000.



### HMGA2 / HMGI-C antibody (pAb) tested by Western blot.

Detection of HMGA2 / HMGI-C by Western blot. Nuclear extract (25 µg) of mouse ES cells probed with HMGI-C / HMGA2 antibody at a dilution of 1:3,000.