

HMGA1 antibody (pAb)

Catalog Nos: 39615, 39650

Isotype: Serum

Application(s): ICC, IF, WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: None

Host: Rabbit

Molecular Weight: 20 kDa

Background: High mobility group (HMG) proteins are nuclear proteins that may significantly affect DNA interactions by altering double-stranded DNA structure. HMGA1 is frequently acetylated and is found in the nucleus.

Immunogen: This HMGA1 antibody was raised against a peptide corresponding to the N-terminus of human HMGA1.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

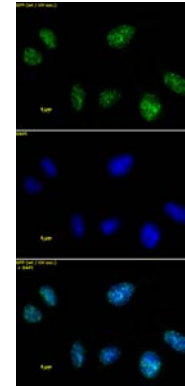
Applications Validated by Active Motif:

ICC/IF: 1:1,000 dilution

WB: 1:500 - 1:2,000 dilution

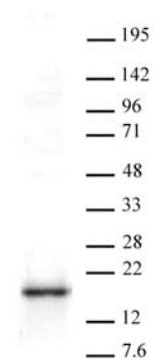
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



HMGA1 antibody tested by immunofluorescence.

Detection of HMGA1 by immunofluorescence. HeLa cells were stained with HMGA1 antibody at a dilution of 1:1,000. Top panel: HMGA1 antibody staining. Middle panel: DAPI. Bottom panel: merge.



HMGA1 antibody tested by Western blot.

HeLa nuclear extract (20 μ g per lane) was probed with HMGA1 pAb (1:500 dilution). Lane 1: no treatment.