

HMGB1 antibody (rAb), 100 µg

Catalog No: 91767**Clone:** 230182G6**Application(s):** WB**Reactivity:** Human, Mouse, Rat**Quantity:** 100 µg**Purification:** Protein A Chromatography**Host:** Rabbit**Isotype:** IgG**Molecular Weight:** 25 kDa

Background: HMGB1, high-mobility group box 1 (also known as high-mobility group protein B1, Amphoterin, HMG-1, HMG1, or SBP-1 belongs to the HMGB family and is involved in unwinding double stranded DNA.

Immunogen: This antibody was raised against the HMGB1 fusion protein.

Buffer: Purified IgG in PBS, pH 7.3, with 50% glycerol and 0.02% sodium azide. Sodium azide is highly toxic.

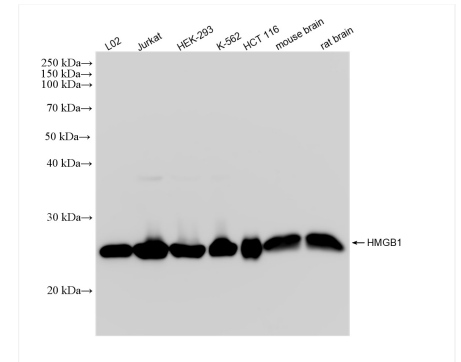
Application Notes:

Applications Validated by Active Motif:

WB: 1:5000-1:50000 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.

**HMGB1 antibody (rAb) tested by Western blot.**

Detection of HMGB1 by Western blot analysis using various lysates and probed with HMGB1 antibody at 1:10000 dilution.