

CBP80 antibody (rAb), 100 µg

Catalog No: 91585**Clone:** 240012G8**Application(s):** WB**Reactivity:** Human**Quantity:** 100 µg**Purification:** Protein A Chromatography**Host:** Rabbit**Isotype:** IgG**Molecular Weight:** 80 kDa

Background: Nuclear cap binding protein subunit 1, 80kDa (NCBP1, synonyms: NCBP, **CBP80**) is a subunit of a cap-binding protein complex (CBC) that consists of a CBP20 subunit, which binds the cap, and a NCBP1 subunit, which ensures high-affinity cap binding. The CBC complex may play a role in pre-mRNA recognition and splicing.

Immunogen: This antibody was raised against the CBP80 fusion protein.

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

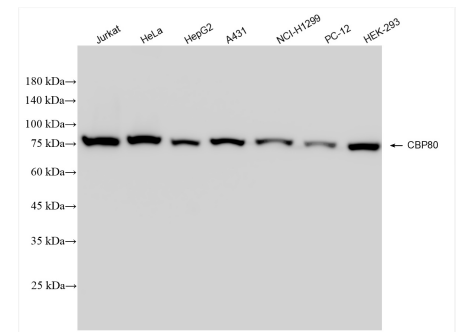
Application Notes:

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.

**CBP80 antibody (rAb) tested by Western blot.**

Detection of CBP80 by Western blot analysis using various lysates and probed with CBP80 antibody at 1:10000 dilution.