

MAX antibody (rAb), 100 µl

Catalog No: 91641**Clone:** 230215000**Application(s):** WB**Reactivity:** Human**Volume:** 100 µl**Purification:** Protein A Chromatography**Host:** Rabbit**Isotype:** IgG**Molecular Weight:** 18 kDa

Background: MAX, a member of MAX family, contains a basic helix-loop-helix (bHLH) domain, is a transcription regulator. MAX can form a sequence-specific DNA-binding protein complex with MYC or MAD which recognizes the core sequence 5'-CAC[GA]TG-3'. The MYC-MAX complex is a transcriptional activator, while the MAD-MAX complex is a repressor. MAX could dimerized with another bHLH protein to form a heterodimer, such MYC or MAD. Thus bands recognized by this antibody much larger than predicted.

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

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

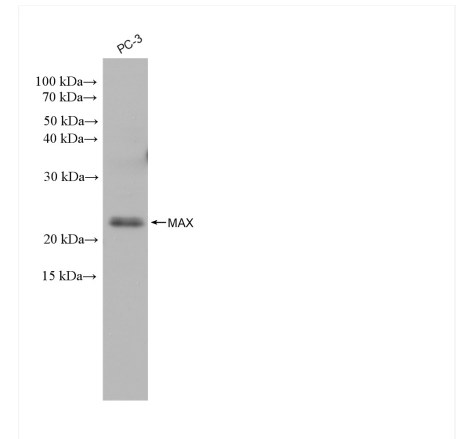
Application Notes:

Applications Validated by Active Motif:

WB: 1:1000-1:8000 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.

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

Detection of MAX by Western blot analysis using 30 µg of PC-3 and probed with MAX antibody at 1:4000 dilution.