

## AIB3 antibody (pAb)

---

**Catalog Nos:** 39835, 39836

**RRID:** AB\_2615018

**Isotype:** IgG

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

**Reactivity:** Human, Mouse

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 250 kDa

---

**Background:** AIB3 (Amplified in Breast Cancer 3, Nuclear Receptor Coactivator 6, NCoA6) is a chromatin-associated protein that interacts with nuclear hormone receptors (antibodies, kits and proteins) to modify their activity as transcriptional activators. AIB3 / NCoA6 was originally isolated as a ligand-dependent nuclear receptor interacting protein. It is a co-activator that directly binds nuclear receptors and stimulates their transcriptional activity in a hormone-dependent fashion. AIB3 is a multifunctional co-activator necessary for transcriptional activation of a wide spectrum of target genes. AIB3 is involved in growth, development, wound healing and maintenance of energy homeostasis. It can enhance the transcriptional activation of a variety of transcription factors and also associates with a number of important regulators of transcription including CBP and p300. AIB3 / NCoA6 is a component of at least three distinct MLL complexes, suggesting that AIB3 plays a fundamental role in transcriptional activation by modulating chromatin structure through histone methylation. AIB3 may also be involved in the co-activation of the NF-kappa-B pathway. It is amplified and overexpressed in breast, colon and lung cancers. AIB3 function is essential in mammalian development as AIB3 null embryos do not survive. Additionally, AIB null spontaneously undergo apoptosis, indicating that AIB3 has anti-apoptotic and pro-survival functions.

**Immunogen:** This AIB3 antibody was raised against a recombinant protein corresponding to amino acids 22-146 of mouse AIB3.

**Buffer:** Purified IgG in 70 mM Tris-HCl (pH 8), 105 mM NaCl, 70 mM glycine, 30% glycerol, and 0.035% sodium azide. Also contains BSA. Sodium azide is highly toxic.

**Application Notes:**

Published Applications:

IHC

WB

See references for more information. Individual optimization may be required.

**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.