

## Atxn7I3 antibody (mAb)

---

**Catalog Nos:** 61339, 61340

**RRID:** AB\_2793597

**Clone:** 2ATX-2B1

**Isotype:** IgG1k

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

**Reactivity:** Human, Mouse

**Quantities:** 100 µg, 10 µg

**Purification:** Protein G Chromatography

**Host:** Mouse

**Concentration:** 1 µg/µl

**Molecular Weight:** 40 kDa

---

**Background:** **Atxn7I3** (ataxin 7-like 3) is a component of the transcription regulatory histone acetylation (HAT) complex SAGA (Spt-Ada-Gcn5-Acetyltransferase), a multiprotein complex that activates transcription by remodeling chromatin and mediating histone acetylation and deubiquitination. Within the SAGA complex, participates in a subcomplex that specifically deubiquitinates both histones H2A and H2B. The SAGA complex is recruited to specific gene promoters by activators such as MYC, where it is required for transcription. Required for nuclear receptor-mediated transactivation. Within the complex, it is required to recruit USP22 and ENY2 into the SAGA complex.

**Immunogen:** This antibody was raised against a synthetic peptide corresponding to amino acid residue(s) 278-296 of human Atxn7I3.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 32 mM glycine, 0.07 mM EDTA, 30% glycerol, and 0.035% sodium azide.

### Application Notes:

Published Applications:

IP

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. Upon receipt, unconjugated antibodies may be stored at -20°C for up to 2 years. Fluorophore- & enzyme-conjugated antibodies should be stored at 4°C. Fluorophore-conjugated antibodies should be protected from light. Keep reagents on ice when not in storage; to avoid repeated freeze/thaw cycles, we recommend aliquoting items that will be stored frozen into single-use fractions prior to freezing. This product is guaranteed for 6 months from date of receipt.

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