

Atxn7 antibody (mAb)

Catalog No: 61347

RRID: AB_2793601

Clone: 3SCA-1C1

Isotype: IgG1k

Application(s): ICC, IF, IHC, WB

Reactivity: Human, Mouse

Quantity: 100 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 100 kDa

Background: **Atxn7** (ataxin-7) acts as component of the STAGA (SPT3-TAF-GCN5 acetylase) transcription coactivator-HAT (histone acetyltransferase) complex. Mediates the interaction of STAGA complex with the CRX (cone-rod homeobox) and is involved in CRX-dependent gene activation.

Immunogen: This antibody was raised against a recombinant protein corresponding to amino acid residue(s) 1-229 of human Atxn7.

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:

ICC/IF

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. 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.