

HIPK2 antibody (mAb)

Catalog Nos: 39677, 39678

RRID: AB_2615043

Clone: 3E9

Isotype: IgG

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

Reactivity: Human

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: 130 kDa

Background: HIPK2 (Homeodomain Interacting Protein Kinase 2) is a serine/threonine kinase present in the nucleus that interacts with homeodomain transcription factors and acts as a co-repressor of transcription. It inhibits cell growth, as well as promotes apoptosis through the phosphorylation and subsequent degradation of the anti-apoptotic factor CTBP1. HIPK2 also phosphorylates CBX4 upon DNA damage and promotes its E3 SUMO-protein ligase activity.

Immunogen: This HIPK2 antibody was raised against a recombinant protein including amino acids 898-1051 of mouse HIPK2.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

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.