

E1A antibody (mAb)

Catalog Nos: 39771, 39772

RRID: AB_2793336

Clone: M73

Isotype: IgG2a

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

Reactivity: Human, Not Species Specific

Quantities: 100 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Background: Adenovirus E1A protein is a transcriptional activator of adenovirus early promoters. This E1A antibody recognizes E1A from cells infected with Adenovirus 2 or Adenovirus 5, and recognizes three major bands from 30 - 50 kDa in molecular weight.

Immunogen: This E1A antibody was raised against a recombinant protein corresponding to full-length E1A from Adenovirus 2.

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.

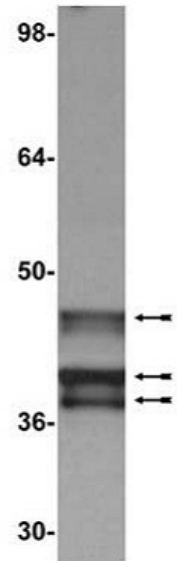
Application Notes:

Applications Validated by Active Motif:

WB: 0.1 - 1 µg/ml 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.



E1A antibody tested by Western blot. HEK293 whole-cell extract probed with E1A antibody (Clone M73) (0.1 µg/ml dilution). Arrows indicate the Ad2 and Ad5 E1A gene products.