

ACTB Monoclonal Antibody

(Catalog # A-0802)

Background

Actin is a highly conserved protein that polymerizes to produce filaments that form cross-linked networks in the cytoplasm of cells. Actin exists in both monomeric (G-actin) and polymeric (F-actin) forms, both forms playing key functions, such as cell motility and contraction. In addition to their role in the cytoplasmic cytoskeleton, G- and F-actin also localize in the nucleus, and regulate gene transcription and motility and repair of damaged DNA.

Description

ACTB Monoclonal Antibody. Unconjugated. Raised in: Mouse

Specificity

Human, Rat, Mouse, Monkey, Dog, Chicken, Hamster, Rabbit, Pig, Sheep

Isotype

IgG

Uniprot ID

P60709

Purification

Affinity Purified

Immunogen

Synthetic Peptide

Formulation

Liquid. PBS, pH 7.4, containing 0.02% sodium azide as Preservative and 50% Glycerol.

Storage

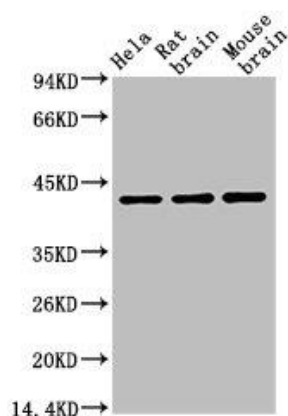
Shipped at 4°C. Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

Cytoplasmic Actin, ACTB, Actin, Actin, Beta, Cytoplasmic 1, B actin, Beta-actin, F actin, β actin, Anti B-actin

Application

ELISA, WB; Recommended dilution: WB:1:500-1:5000



Western blot analysis of HeLa (1), Rat brain (2), Mouse brain (3), (A02) diluted at 1:10000.