

Myc-Tag antibody (mAb)

Catalog Nos: 39279, 39280

RRID: AB_2793216

Clone: 4E12

Isotype: IgG1

Application(s): WB

Reactivity: Mouse, Not Species Specific

Quantities: 200 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

Background: Myc-Tag is a sequence derived from the human c-Myc protein that has been developed into an effective epitope for detection of tagged proteins expressed *in vivo*. Clone 4E12 exhibits superior detection of Myc-Tagged proteins compared to the more common 9E10 clone.

Immunogen: This Myc-Tag antibody was raised against peptide corresponding to amino acids 410-419 of human c-Myc (EQKLISEEDL).

Buffer: Purified rabbit IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

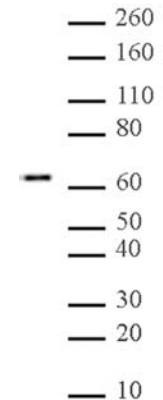
WB: 0.5 - 2 µg/ml dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

This antibody is also available as an AbFlex® engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91203.

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.



Myc-Tag mAb tested by Western blot.

Whole-cell extract (20 µg) of NIH/3T3 cells expressing a Myc-Tagged PP2A-alpha subunit blotted with Myc-Tag mAb (0.5 µg/ml).