

# [K(Ac)40] $\alpha$ -Tubulin polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 6

[View Online »](#)

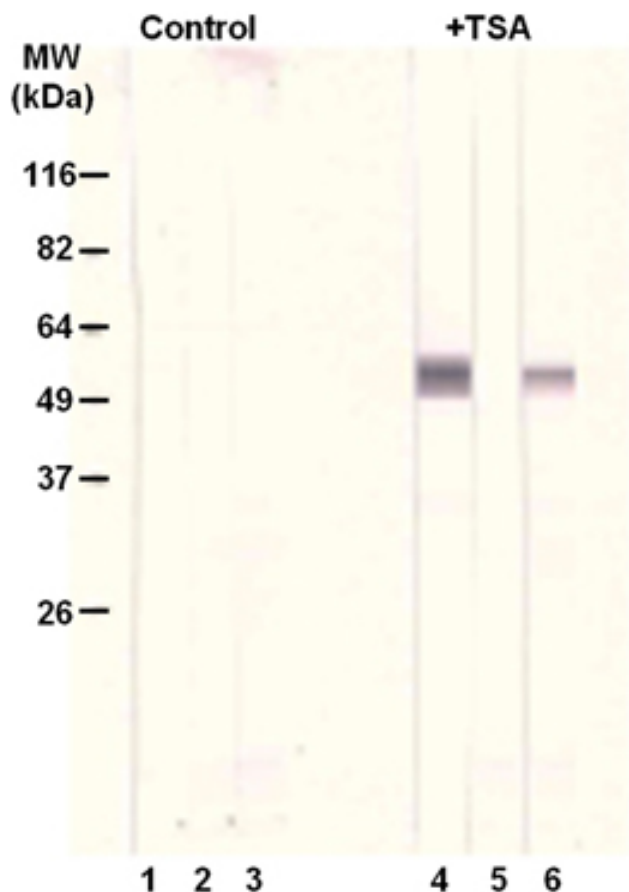
## Ordering Information

[Order Online »](#)

BML-SA452-0100	100 $\mu$ l
----------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Treatment of A549 cells with the histone deacetylase inhibitor trichostatin A (TSA, 0.3  $\mu$ M, 7 hr.) induces acetylation of alpha-tubulin-Lys40. Western blot analysis was performed with A549 cell extracts (100,000 cells per lane) from either vehicle treated (Control) or TSA-treated (+TSA) cells. All lanes were probed with a 1:1000 dilution of the primary antibody, anti-alpha-tubulin (K(Ac)40) (Prod. No. BML-SA452). For lanes 2 and 5 the primary antibody was preincubated with the immunogen peptide (Ac-QMPSPD-K(Ac)-TIGGG[CNH<sub>2</sub>]) and for lanes 3 and 6 it was preincubated with non-acetylated peptide Ac-QMPSPDKTIGGG[C-NH<sub>2</sub>] (peptides: 0.5  $\mu$ g/ $\mu$ l, antibody: 1:100 dilution, 37°C, 1hr.). Secondary antibody for all lanes was donkey anti-rabbit-AP (1:2000; Jackson ImmunoResearch) and the nitrocellulose membrane was developed with BCIP/NBT.

## Handling & Storage

**Handling** Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** WB

**Application Notes** Detects a band of ~50kDa by Western blot.

**Formulation** Liquid. In PBS.

**Host** Rabbit

**Immunogen** Synthetic acetylated peptide (K(Ac)40) corresponding to aa 35-45 (Q35MPSD-K(Ac)-TIGGG45) of human  $\alpha$ -tubulin.

**Recommendation Dilutions/Conditions** Western Blot (1:1,000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Source** Purified from rabbit serum.

**Species Reactivity** Chicken, Primate, Rodent, Xenopus, Zebrafish

**UniProt ID** Q71U36 (TUBA1A), P68363 (TUBA1B), P0DPH7 (TUBA3C)

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#)

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)