

Acetylated α -Tubulin (K40) Recombinant monoclonal antibody (RM318)

The acetylated residue of α -tubulin is K40, which is catalyzed by α -tubulin acetyl-transferase (α -TAT) in humans. The acetylation of K40 on alpha tubulin is a hallmark of stable microtubules.

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-ABS676-0100	100 μ l
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ICC, IF, WB

Clone RM318

Formulation Liquid. In 1x PBS containing 50% glycerol, 1% BSA and 0.09% sodium azide

Host Rabbit

Immunogen Synthetic acetyl-peptide corresponding to the K40 region of human alpha tubulin.

Isotype IgG

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot: 1:1,000-1:2,000IF/ICC: 1:1,000-1:5,000Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from animal origin-free supernatant.

Species Reactivity Human

UniProt ID P68366

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



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-
eu@enzolifesciences.com](mailto:info-eu@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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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