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

100µl

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/ConditionsWestern 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



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com