

# TPPP monoclonal antibody (17C6)

TPPP plays a role in the polymerization of tubulin into microtubules, microtubule bundling and the stabilization of existing microtubules, thus maintaining the integrity of the microtubule network. Plays a role in mitotic spindle assembly and nuclear envelope breakdown.

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

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 1 year after receipt when stored at -20°C.
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Tubulin polymerization-promoting protein, p25
<b>Application</b>	ICC, IHC (PS), WB
<b>Clone</b>	17C6
<b>Formulation</b>	Liquid. In PBS containing 10% glycerol and 0.02% sodium azide.
<b>Host</b>	Rat
<b>Immunogen</b>	Full length recombinant TPPP.
<b>Isotype</b>	IgG2a
<b>Purity Detail</b>	Protein G affinity purified.
<b>Recommendation Dilutions/Conditions</b>	Immunocytochemistry (10µg/ml)Immunohistochemistry (paraffin sections, 10µg/ml)Western Blot (200ng/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Source</b>	Purified from concentrated hybridoma tissue culture supernatant.
<b>Species Reactivity</b>	Human, Monkey, Mouse

**Specificity**

Recognizes full-length endogenous TPPP/p25.

**UniProt ID**

O94811

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

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