

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 »](#)

Ordering Information

[Order Online »](#)

ALX-803-342-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Tubulin polymerization-promoting protein, p25
Application	ICC, IHC (PS), WB
Clone	17C6
Formulation	Liquid. In PBS containing 10% glycerol and 0.02% sodium azide.
Host	Rat
Immunogen	Full lenght recombinant TPPP.
Isotype	IgG2a
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	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.
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	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