
TPPP Monoclonal Antibody [17C6]

(Catalog No. A-0520)

Background

Tubulin polymerization-promoting protein (TPPP) is a brain-specific protein expressed mainly in myelinating oligodendrocytes of the central nervous system. It is thought to play a role in the polymerization of tubulin into microtubules, microtubule bundling and the maintenance of existing microtubules, thereby conserving the integrity of the microtubule network. TPPP may also be involved in mitotic spindle assembly and nuclear envelope breakdown.

Description

Rat monoclonal antibody to TPPP, Clone 17C6

Formulation

Liquid. In PBS containing 10% glycerol and 0.02% sodium azide

Immunogen

Full length recombinant TPPP

Specificity

Human, Mouse, Monkey

Isotype

IgG2a

Purification

Protein G-affinity purified

Storage

Store at +4°C for short term. Aliquot and store at -20°C for long term. Avoid multiple freeze/thaw cycles.

Application

ICC (10 µg/ml), IHC (paraffin sections 10 µg/ml), WB (200 ng/ml)

Ordering Information

Products	Size	Cat. No.
TPPP Monoclonal Antibody [17C6]	100 µg	A-0520-100

This product is for research purposes only. Not intended for use in diagnostic procedures.