β-Tubulin Recombinant monoclonal antibody (S11B) (Rabbit IgGλ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS614-0200

200µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.

Handling Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name TUBB

Application ELISA, WB

Clone S11B

Crossreactivity No crossreactivity with alpha-tubulin.

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Rabbit

Immunogen Human MBP (microtubule-binding protein).

lsotype lgGλ

Purity Detail Protein A affinity purified.

Species Reactivity Human

Technical Info / Product NotesThis full-length, chimeric rabbit antibody was made using

the variable domain sequences of the original Human scFv format, for improved compatibility with existing reagents,

assays and techniques.

UniProt ID Q13509

Last modified: May 29, 2024



info-

eu@enzolifesciences.com