

β III-Tubulin (human) monoclonal antibody (TU-20)

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

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Class III β -tubulin
Application	ELISA, Flow Cytometry, ICC, IHC (FS), IHC (PS), WB
Clone	TU-20
Formulation	Liquid. In PBS, pH 7.4, containing 15 mM sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to aa 441-448 (E441SESQGPK448) of human β III-tubulin (class III β -tubulin).
Isotype	IgG1
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections, 10 μ g/ml; pretreatment with 0.1% pepsin (trypsin) in 0.1M HCl; incubation for 30 min. at room temperature or high temperature citrate buffer antigen retrieval)Western Blot (1-2 μ g/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Specificity	Recognizes β III-tubulin located in neurons.

Technical Info / Product Notes

Application Note:

[Image-Based Analysis of a Human Neurosphere Stem Cell Model for the Evaluation of Potential Neurotoxicants](#)

UniProt ID

Q13509

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)