

Tau (human) monoclonal antibody (Tau 12)

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

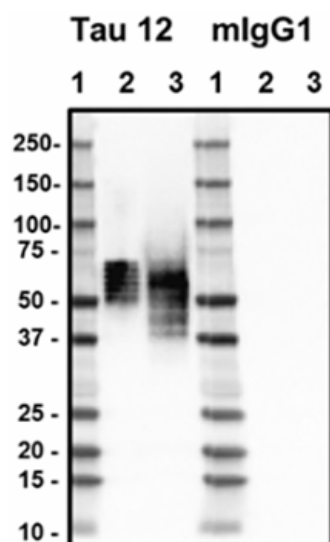
Ordering Information

[Order Online »](#)

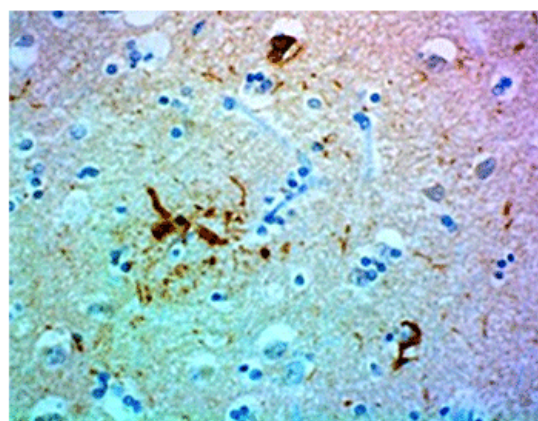
| | |
|-----------------|-------|
| ENZ-ABS216-0100 | 0.1ml |
| ENZ-ABS216-0500 | 0.5ml |

Manuals, SDS & CofA

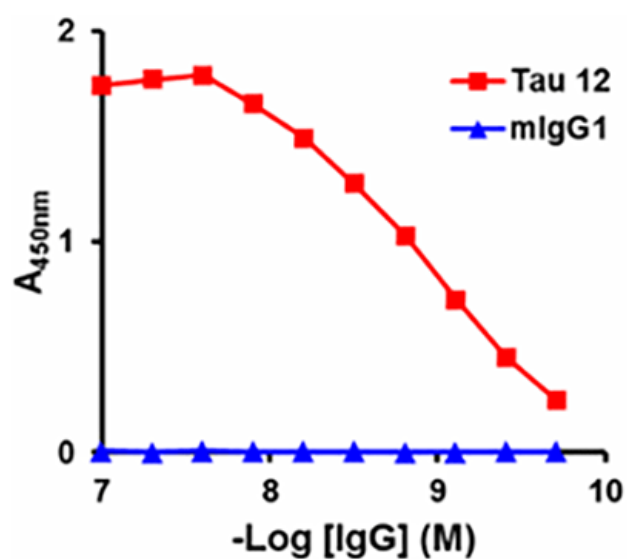
[View Online »](#)



Western blot of Tau (human) monoclonal antibody (Tau 12). Lane 1: Molecular weight marker; Lane 2: Tau protein ladder; Lane 3: 20 µg of human brain lysate.



IHC staining of Tau (human) monoclonal antibody (Tau 12) on formalin-fixed paraffin-embedded human Alzheimer's disease brain tissue.



Direct ELISA of Tau (human) monoclonal antibody (Tau 12) and isotype-matched control mouse IgG1 binding to plate-immobilized recombinant human 2N4R Tau protein.

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

| | |
|----------------------|---|
| Alternative Name | Microtubule-associated protein Tau |
| Application | ELISA, IHC, WB |
| Clone | Tau 12 |
| Formulation | Liquid. In PBS. Does not contain any preservatives or carrier proteins. |
| Host | Mouse |
| Isotype | IgG1 |
| Purity Detail | Purified. |
| Species Reactivity | Human |
| Specificity | Recognizes an epitope between aa 6-18 of human Tau. |
| UniProt ID | P10636 |
| 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-
fr@enzolifesciences.com](mailto:info-fr@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)