

Tau-383 (0N4R variant) (human), (recombinant)

The tau proteins are major neuronal microtubule-associated proteins that function in the stabilization of microtubules and accumulate in various abnormal filaments including neurofibrillary tangles, neuropil threads and amyloid plaques in the pathology of Alzheimer's disease. Six tau isoforms are derived from alternative mRNA splice variants that originate from a single gene and result in mature proteins that vary in size from 352 to 441 amino acids (36.8 to 45.9 kDa). The six isoforms differ from one another in having three or four microtubule binding repeats (R) and two, one or zero amino terminal inserts (N). While the fetal brain contains a single isoform of tau (Tau-352), the adult brain has several isoforms.

Ordering Information

[Order Online »](#)

BML-SE324-0100	100µg
-----------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

This protein is purified without any acid treatment and is suitable for promoting microtubule assembly, and for hyperphosphorylation-induced self-assembly into filaments.

Formulation

Lyophilized from 50mM MES, pH 6.8, 100mM NaCl and 0.5mM EGTA.

MW

40.0 kDa

Source

Produced in *E. coli*.

UniProt ID

P10636

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)