

Tau dGAE (297-391) Monomers

Human Recombinant Tau dGAE (AA297-391)
Monomers
Catalog No. SPR-501



Discovery through Partnership | Excellence through Quality

EAGLE BIOSCIENCES, INC.
20A NW BLVD, SUITE 112 NASHUA, NH 03063
P: 617-419-2019 F: 617-419-1110
WWW.EAGLEBIO.COM — INFO@EAGLEBIO.COM



Product Name

Tau dGAE (297-391) Monomers

Description

Human Recombinant Tau dGAE (AA297-391) Monomers

Applications

WB, SDS PAGE, In vitro Assay

Concentration

Lot/batch specific. See included datasheet.

Conjugates

No tag

Nature

Recombinant

Species

Human

Expression System

E. coli

Amino Acid Sequence

IKHVPGGGSVQIVYKPVLDLSKVTSKCGSLGNIHHKPGGGQVEVKSEKLDKDRVQSKIGSLDNITHVPGGGNKKIETHKLTFRN
AKAKTDHGAE

Purity

>95%

Other Resources

Protein Length

Fragment of full length wild-type Tau 2N4R (297 - 391aa)

Field Of Use

Not for use in humans. Not for use in diagnostics or therapeutics. For in vitro research use only.

Properties

Storage Buffer

10mM PB pH 7.4, 10mM DTT

Storage Temperature

-80°C

Shipping Temperature

Dry Ice. Shipping note: Product will be shipped separately from other products purchased in the same order.

Purification

Ion-exchange, ammonium sulfate precipitation and SEC purified

Cite This Product

Human Recombinant Tau dGAE (297-391) Monomers (StressMarq Biosciences Inc., Victoria BC CANADA, Catalog # SPR-501)

Certificate Of Analysis

Protein certified >95% pure on SDS-PAGE & Nanodrop analysis. Low endotoxin <5 EU/mL @ 2mg/mL.

Biological Description

Alternative Names

Tau monomer, Tau protein monomer, Tau protein, microtubule-associated protein Tau, MAPT, MAP, microtubule-associated protein, Truncated Tau Protein Monomer, Paired Helical Filament-Tau, Phf-Tau, Neurofibrillary Tangle Protein, Tau dGAE Protein, dGAE

Research Areas

Alzheimer's Disease, Axon Markers, Cell Markers, Cell Signaling, Cytoskeleton, Microtubules, MT Associated Proteins, Neurodegeneration, Neuron Markers, Neuroscience, Tangles & Tau

Swiss Prot

P10636-8

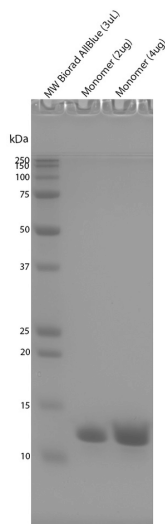
Scientific Background

Filamentous tau inclusions are a hallmark of many neurodegenerative diseases, including Alzheimer's disease (AD) and Chronic Traumatic Encephalopathy (CTE), collectively called tauopathies. Advances in Cryo-EM have revealed that tau filaments isolated from individuals with a particular neurodegenerative disease share a distinct tau fold – i.e. an AD-isolated Tau filaments' fold is distinct from a CTE-isolated Tau filaments' fold (1-3). Utilizing Tau filaments with the correct disease-specific fold is an important goal towards better mimicking specific human diseases in cellular and in vivo models. Recent Cryo-EM studies have demonstrated that recombinantly generated Tau dGAE monomers will form the disease-isolated AD or CTE Tau filament folds under highly specific conditions in vitro (4, 5). StressMarq's SPR-501 Tau (297-391) dGAE monomers are purified following these exact published procedures and can be utilized to form these distinct folds using specific aggregation conditions.

References

1. Goedert, Eisenberg and Crowther. 2017. Propagation of Tau Aggregates and Neurodegeneration. *Annu Rev Neurosci*. DOI: <https://doi.org/10.1146/annurev-neuro-072116-031153>
2. Fitzpatrick et al. 2017. Cryo-EM structures of tau filaments from Alzheimer's disease. *Nature*. DOI: [10.1038/nature23002](https://doi.org/10.1038/nature23002)
3. Falcon et al. 2019. Novel tau filament fold in chronic traumatic encephalopathy encloses hydrophobic molecules. *Nature*. DOI: <https://doi.org/10.1038/s41586-019-1026-5>.
4. Lovestam et al. 2022. Assembly of Recombinant Tau into Filaments Identical to those of Alzheimer's disease and Chronic Traumatic Encephalopathy. *eLife*. DOI: <https://doi.org/10.7554/eLife.76494>
5. Lovestam et al. 2023. Disease-specific Tau Filaments Assemble via Polymorphic Intermediates. *bioRxiv*. <https://doi.org/10.1101/2023.07.24.550295>

Product Images



SDS-PAGE of purified E. coli expressed hTau (dGAE) (SPR-501) on a 12% Tris-Glycine gel.



TEM of Tau dGAE AD-mimic fibrils (SPR-502) generated from Tau dGAE monomers (SPR-501) by shaking 200 rpm at 37°C for 48 hours in 10 mM PB 10 mM DTT pH 7.4 with 200 mM MgCl₂ added (Lovestam et al. 2022, eLife). Negative stain transmission electron microscopy images acquired at 80 Kv on carbon coated 400 mesh copper grids using phosphotungstic acid and uranyl acetate stain. Scale bar = 200 nm.

Product Citations (0)

Currently there are no citations for this product.

Reviews

There are no reviews yet.