

# Tau-383 (0N4R) Wild-Type Monomers

Human Recombinant Tau-383 (0N4R) Wild-Type  
Monomers  
Catalog No. SPR-509



Discovery through Partnership | Excellence through Quality

distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com) info@eaglebio.com



## Product Name

Tau-383 (0N4R) Wild-Type Monomers

## Description

Human Recombinant Tau-383 (0N4R) Wild-Type 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

MAEPRQEFVEMEDHAGTYGLGDRKDQGGYTMHQDQEGD TDAGLKAE EAGIGDTPSLEDEAAGHVTQARMVSKSKDGTGS  
DDKKAKGADGKTKIATPRGAAPPGQKGQANATRIPAKTPPAPKTPPSSGEPKSGDRSGYSSPGSPGTPGSRSRTPSLTPPTR  
EPKKVAVVRTPPKSPSSAKSRLQTAPVPMPDLKNVSKIGSTENLKHQPGGKQVQIINKKLDLSNVQSKCGSKDNIKHVPGGGS  
VQIVYKPV DLSKVT SKCGSLGNIHHKPGGGQVEVKSEKLDKDRVQSKIGSLDNITHVPGGGNKKIETHKLTFR ENAKAKTDHG  
AEIVYKSPVWSGDTPRHLSNVSS TGSIDMV DSPQLATLADEV SASLAKQGL

## Purity

---

>95%

---

## Other Resources

---

## Protein Length

---

383 aa

---

## Protein Size

---

40.007 kDa

---

## 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 Hepes pH 7.4, 100mM NaCl

---

### 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 Purified

---

### Cite This Product

---

Human Recombinant Tau-383 (0N4R) Wild-Type Monomers (StressMarq Biosciences Inc., Victoria BC CANADA, Catalog # SPR-509)

---

### Certificate Of Analysis

---

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

---

### Other Relevant Information

---

For corresponding PFFs, see catalog# SPR-510

---

## Biological Description

---

### Alternative Names

Tau-D, Tau 383

---

### Research Areas

Alzheimer's Disease, Neurodegeneration, Neuroscience, Tangles & Tau

---

### Swiss Prot

P10636-6

---

### Scientific Background

Several tau isoforms, including 0N4R, are expressed in the human brain, and the existence of multiple human tauopathies with distinct fibril morphologies suggests different molecular conformers incorporating different isoforms may exist (1, 2). NMR data indicates that both 3R and 4R tau are incorporated into AD-tau seeded fibrils (3).

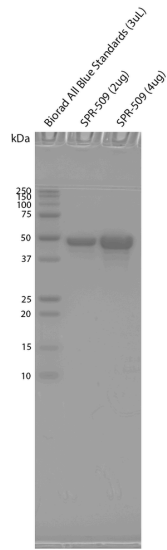
---

### References

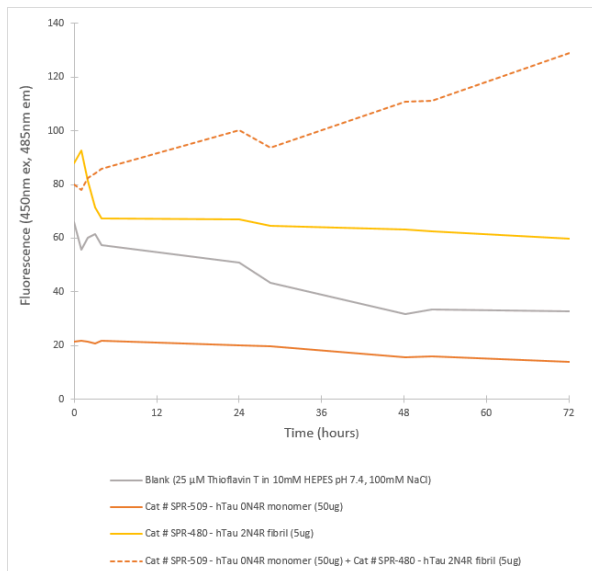
1. Goedert et al. 1989. Multiple isoforms of human microtubule-associated protein tau: sequences and localization in neurofibrillary tangles of Alzheimer's disease. *Neuron*. DOI: 10.1016/0896-6273(89)90210-9
  2. Goedert, Eisenberg and Crowther. 2017. Propagation of Tau Aggregates and Neurodegeneration. *Annual Review of Neuroscience*. DOI: 10.1146/annurev-neuro-072116-031153
  3. Dregni et al. 2022. Fluent molecular mixing of Tau isoforms in Alzheimer's disease neurofibrillary tangles. *Nature communications*. DOI: 10.1038/s41467-022-30585-0
- 

## Product Images

---



SDS-PAGE analysis of purified E. coli expressed human tau-383 (0N4R) monomers on a 12% Tris-Glycine gel.



In vitro seeding activity of human tau-383 (0N4R) monomers in ThT assay. Full length, wild-type tau 2N4R fibrils (SPR-480) seed fibril formation of human Tau-383 (0N4R) monomers over 72 hours. Reactions (100uL) shaken at 600 rpm in Greiner-Bio 96 Well Non-Binding Cell Culture Microplates, Black (Greiner-Bio Catalog #655900) at 37oC and read with an XPS Microplate Reader set at 450nmex/485nmem.

## Product Citations

## Reviews

---

There are no reviews yet.