



Product Data Sheet

Product Name:	Recombinant Human Tau (Tau-441), Biotin Labeled
Catalog Number:	AS-56084
Lot Number:	See label on vial
Amount:	50 µg
Source:	The recombinant human Tau (Tau-441) (GenBank Accession # AAC04279.1) was expressed and purified from <i>E. coli</i> and conjugated with biotin.
Purity:	Greater than 90% as determined by SDS-PAGE.
DOS:	See label on the vial
Storage:	Biotinylated human Tau (Tau-441) is supplied frozen at 1 mg/ml in 10 mM sodium phosphate buffer (pH=7.4). Store at 2-4 °C for immediate use within 1 week or at -80 °C for long-term storage. Avoid repeated freeze-thaw cycles.
Description:	<p>Microtubule associated protein (Tau) is found predominantly in the central neural system and its major function is to promote assembly and to stabilize neuronal microtubules.¹⁻⁵ Six isoforms of Tau were identified in humans that are differentiated by the exclusion or inclusion of exons 2, 3, and 10.²⁻⁵ Tau-441 is the longest of Tau isoforms that consists of 441 amino acids with molecular mass of 45.8 kDa.</p> <p>Tau441-Biotin conjugates have been reported in applications such as Tau species identification⁶ and different binding assays⁷⁻¹⁰.</p>

Related Products

Product Name	Cat. #
Recombinant Human Tau (Tau-441) Protein	AS-55556
Recombinant Human Tau (Tau-441) Protein, GST tagged	AS-55557

References:

1. Bulic B., Neuropharmacology. 59 (2010):276-89.
2. Voss K., Mol. Neurodegen. 4 (2009):1-12.
3. Rankin CA., Mol. Neurodegen. 2 (2007):1-14.
4. Patterson KR., J. of Biol. Chem. 286 (2011):23063-76.
5. Wang Y., Biochem. Soc. Trans. 38 (2010):955-61.
6. Derisbourg M., Sci Rep. 5 (2015):9659.
7. Stopschinski BE., J Biol Chem. 295 (2020):2974-83.
8. Berrocal M., Biochim Biophys Acta. 1852 (2015):1465-76.
9. Flach K., Biochim Biophys Acta. 1842 (2014):1527-38.
10. Makrides V., J Biol Chem. 278 (2003):33298-304.

For in vitro research use only.