

## Product Information

<b>Description:</b>	Recombinant human full-length Tau-441, also known as Tau-F. This construct does not contain a tag but includes the mutation of interest P301L.
<b>Species:</b>	Human
<b>Construct:</b>	Tau-441 (P301L) (Full Length)
<b>Mutation:</b>	P301L
<b>Concentration:</b>	0.20 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 0.25 mM DTT, 0.1 mM PMSF, and 25% glycerol
<b>MW:</b>	64 kDa
<b>Uniprot:</b>	P10636-8
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining

