



## Product Data Sheet

---

Product Name:	<b>Anti-Tau (paired262) Blocking Peptide</b>
Catalog Number:	54974P
Lot Number:	See label on vial
Product Description:	Antibody blocking peptide supplied with the same amino acid sequence as the immunizing peptide used for the antibody production of anti-TAU (paired262), cat# 54974. Each package contains 100 µg peptide conjugated to BSA in 100 µl of 1xPBS (1 mg/ml).
Quality Control:	The quality of the peptide was evaluated by HPLC and mass spectrometry.
Application Notes:	Blocking peptide may be used as a control in applications such as ELISA, western blot (WB), immunohistochemistry (IHC), immunofluorescence (IF), flow cytometry (FC), immunocytochemistry (ICC) and immunoprecipitation (IP) by blocking the binding of specific antibody to the target antigen. For blocking purpose, the antibody can be incubated with the blocking peptide at optimal concentrations for 2 hours at room temperature or overnight at 4 °C in selected buffer (generally the buffer used for primary antibody incubation). The ratio of the blocking peptide and antibody can be 5:1 to 10:1 or higher (w/w).
Storage:	Store at -20 °C for up to 12 months upon receiving the product.

This product is for *in vitro* research use only.