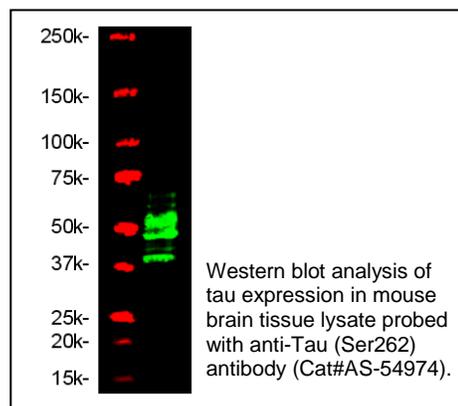


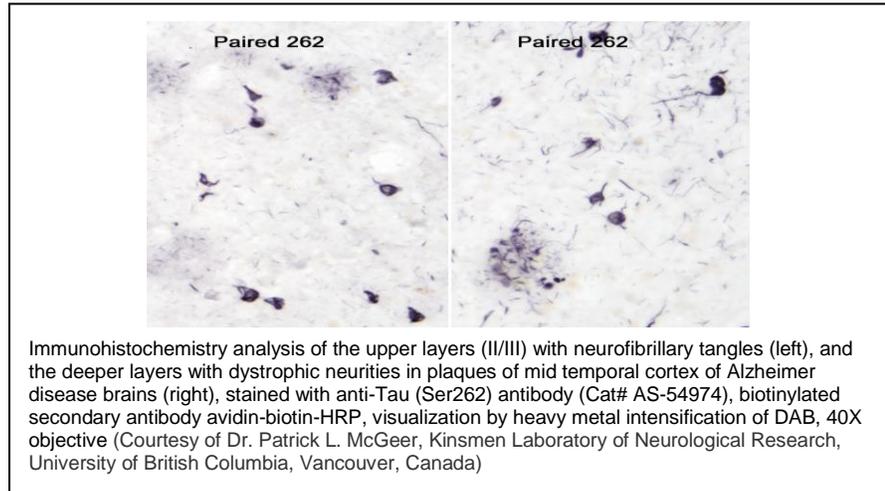
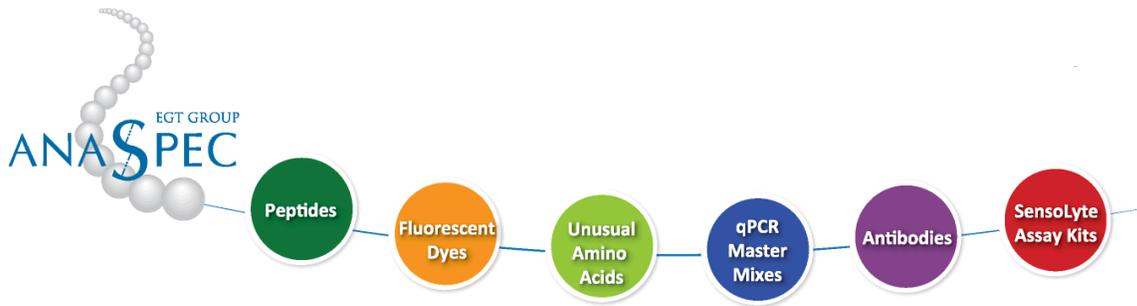
Update: July, 2017

Product Data Sheet

Product Name:	Anti-Tau (Ser262) <i>Rabbit Polyclonal Antibody</i>
Catalog Number:	AS-54974
Lot Number:	See label on vial
Storage Buffer:	1X PBS (pH 7.4) containing 0.05% sodium azide < 0.1% BSA
Size:	50 µg
Concentration:	0.2 mg/mL
Immunogen:	KLH conjugated with synthetic peptides corresponding to human Tau surrounding Serine 262 region
Species Reactivity:	Species reactivity includes human, mouse, rat and bovine while others remain unknown. The antibody was evaluated for specificity by dot blot, Western blot and immunohistochemistry.
Application Notes:	The following concentration ranges are recommended starting points for this product. Optimal working concentrations should be determined by the investigator for specific applications.

Dot blot:	0.5-2.0 µg/mL
Western blot:	0.5-2.0 µg/mL
Immunohistochemistry:	5.0-20.0 µg/mL





Background:

Tau is a collection of microtubule-associated proteins that is involved in microtubule assembly and stabilization.¹ In adult human brain, 6 isoforms ranging between 352 and 441 amino acids in length are produced as a result of alternative RNA splicing.^{2,3} The expression of tau isoforms is developmentally regulated, as only the smallest tau polypeptide is expressed in fetal brain. Hyperphosphorylated Tau is the major component of the paired helical filament of Alzheimer's disease. Phosphorylation-dependent anti-Tau antibodies are used to identify specific amino acids that are phosphorylated in tau from normal brain and Alzheimer's disease brain. The Tau proteins, esp. in developing brain and in Alzheimer brains, were found to phosphorylate *in vivo* at many different sites.⁴

Storage:

Store at 4°C for 1-2 weeks. Aliquot and store at -20°C up to 1 year. Avoid freeze and thaw cycle.

References:

1. Cleveland DW, et al (1977) *J. Mol. Biol.* **116**, 207-225
2. Goedert M, et al (1989) *Neuron* **3**, 519-526.
3. Geodert M, et al (1989) *EMBO J.* **8**, 393-399.
4. Billingsley M et al (1997) *Biochem J* **323**, 577-591.

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