



Tau [pS⁴⁰⁰] Phospho and Non-Phospho Peptide Sample Pair

PRODUCT ANALYSIS SHEET

Catalog Number:	04-754Z		
Lot Number:			
Components:	Part Number	Quantity	Lot Number
	04-754P	0.1 mg	
	04-754N	0.1 mg	
Description:	Positive and negative control peptides for the Tau phosphoserine ⁴⁰⁰ rabbit polyclonal antibody (Cat. # 44-754)		
Data for phospho peptide:	04-754P		
Molecular Weight	1674		
Purity	>90% by HPLC analysis (purity based on peak area)		
Peptide Content	80 ± 4%		
Data for non-phospho peptide:	04-754N		
Molecular Weight	1594		
Purity	>85% by HPLC analysis (purity based on peak area)		
Peptide Content	80 ± 4%		
Supplied As:	Trifluoroacetate salt; 0.1 mg peptide/vial		
Physical Appearance:	Lyophilized powder		
Solubility/Reconstitution:	We recommend that each vial be briefly centrifuged prior to opening to bring the contents to the bottom. To reconstitute the peptides, pipette 100 µL sterile, distilled H ₂ O into each vial. The peptides may be gently vortexed or triturated at room temperature to ensure complete solubilization. If the peptides are found to be slow to dissolve, the pH of the solutions may be neutralized by addition of small volumes (1-2 µL) 1.0 N ammonium hydroxide; alternatively, the peptides may be solubilized by addition of small volumes of organic solvent, such as ethanol or DMSO.		
Storage:	-20°C		
Expiration Date:	Expires one year from date of receipt when stored as instructed.		

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI 04754Z

(Rev 08/09) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.