

PRODUCT DATA SHEET

Revised: 18 April 2005

Printed: 29 January 2009; page 1 of 1



Product name(s):	Rpn11 blocking peptide for rabbit pAb PW9625
-------------------------	---

Catalogue number:	PP9630	Batch number:	XXXX	Expiry date:	12 months from receipt
--------------------------	--------	----------------------	------	---------------------	------------------------

Product information:

Rpn11, a non-ATPase subunit situated in the lid subcomplex of the 19S regulator, partakes in the processing of substrates tagged with polyubiquitin chains or in their removal from substrates bound to the proteasome. Rpn11 also plays a role in maintaining mitochondrial integrity, tubular structure and proper function. The recent finding that Rpn11 participates in proteasome-associated deubiquitylation focuses interest on the MPN+ (Mpr1, Pad1, N-terminal)/JAMM (JAB1/MPN/Mov34) metalloprotease site in its N-terminal domain¹.

The peptide corresponding to amino acid residues 19-32 of Rpn11 (accession number NP_005796), was synthesised *via* a manual Fmoc/t-butyl-based solid-phase synthesis strategy. Purification was carried out by preparative high performance liquid chromatography (hplc). The purified product was analysed by reverse-phase hplc, by amino acid analysis following acid hydrolysis, and by mass spectrometry. Peptide PP9630 is supplied for control purposes for use with antibody PW9625.

Analytical and physico-chemical data:

Amino acid analysis:	Not determined.
Purity:	Determined to be $\geq 95\%$ by hplc.
Spectroscopic analysis:	Laser desorption mass spectrometry shows molecular ion at 1579.6 Da being consistent with the desired product.
Form:	Lyophilised solid supplied as gross peptide.
Solubility:	Soluble in aqueous solution.

Stability, storage and specific hazard data:

Store solid cold and dry at -20°C . Store solutions at -20°C for up to three months.

References:

1. Rinaldi, T., Elah, P., Gambadoro, A., Zilli, S., Maytal-kivity, V., Frontali, L. and Glickman M. Participation of the proteasomal lid subunit Rpn11 in mitochondrial morphology and function is mapped to a distinct C-terminal domain. *Biochem. J.*, **381**, 275–285 (2004).