

PRODUCT DATA SHEET

Revised: 23 March 2006
Printed: 28 January 2009; page 1 of 1



Product name(s):	Bz-Val-Gly-Arg-AMC				
-------------------------	---------------------------	--	--	--	--

Catalogue number:	BW 9375	Batch number:	Z05953	Expiry date:	12 months from receipt
--------------------------	---------	----------------------	--------	---------------------	------------------------

Product information:
A substrate for the trypsin-like activity associated with the 20S proteasome. Excitation: 380nm, emission: 460 nm.

Structure:
Bz-Val-Gly-Arg-AMC.TFA Bz = benzoyl; AMC = 7-amido-4-methylcoumarin; TFA = trifluoroacetate salt

Molecular formula:	$C_{30}H_{37}N_7O_6.TFA$
Molecular weight:	591.67Da (excluding trifluoro-acetate salt)
CAS number:	Not allocated

Analytical and physico-chemical data:	
Melting point:	138°C
Purity:	>99% as judged by high performance and thin layer chromatography.
Spectroscopic analysis:	Mass spectrum is fully consistent with anticipated molecular weight.
Form:	White powder.
Solubility:	Soluble in methanol or DMSO.

Stability, storage and specific hazard data:
Store solid cold and dry at -20C; stable for at least one year. Store solutions at -20C for up to 1 month.

References:
1. Andersson, M., Sjostrand, J., Karlsson, J.O. Differential inhibition of three peptidase activities of the proteasome in human lens epithelium by heat and oxidation. <i>Exp Eye Res.</i> , 69 (1):129-138 (1999)