

Product Information Sheet

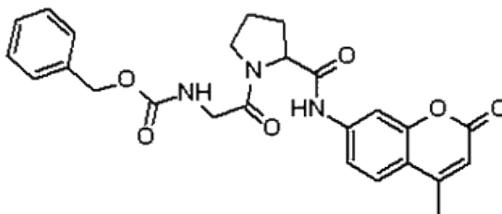
Ordering Information

Product Number: 13477
Product Name: Z-Gly-Pro-AMC
Unit Size: 5 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 463.49

Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 351 nm
Emission Wavelength: 430 nm

Application Notes

Z-Gly-Pro-AMC is a fluorogenic substrate for the determination of post-proline cleaving enzyme (prolyl endopeptidase). This substrate is hydrolyzed by prolyl endopeptidase to generate highly fluorescent 7-amido-4-methylcoumarin (AMC).