

Product Information Sheet

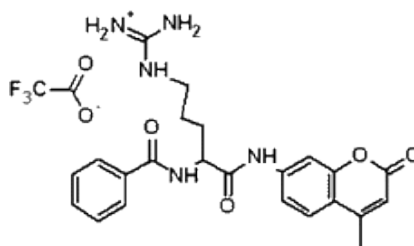
Ordering Information

Product Number: 13478
Product Name: Benzoyl-Arg-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: 549.50

Chemical Structure:



Soluble in: DMSO

Excitation Wavelength: 351 nm

Emission Wavelength: 430 nm

Application Notes

This Benzoyl-Arg-AMC is a fluorogenic aminopeptidase and endopeptidase substrate used for the detection of enzymatic activity of human alpha-thrombin. The non-fluorescent substrate generates a bright blue fluorescent AMC product that has Ex/Em = 351/430 nm, and can be easily detected with the DAPI filter set.