

Catalog number: 13461
Unit size: 5 mg

Component	Storage	Amount (Cat No. 13461)
BOC-Val-Pro-Arg-AMC	Freeze (< -15 °C), Minimize light exposure	1 vial (5 mg)

KEY PARAMETERS

Instrument: Fluorescence microplate reader
Excitation: 360 nm
Emission: 470 nm
Recommended plate: Solid black

PREPARATION OF STOCK SOLUTIONS

1. BOC-Val-Pro-Arg-AMC stock solution

Prepare a 5 to 10 mM stock solution in DMSO. e.g. To prepare 10 mM stock solution, mix 1 mg of powder to 135 μ L DMSO.

Note: Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles

PREPARATION OF WORKING SOLUTION

Note: Prepare the BOC-Val-Pro-Arg-AMC working solution fresh before each experiment, and protect it from light.

Component	Volume
Substrate stock solution (10 mM)	25 µL
DTT (1M)	100 µL
EDTA (100 mM)	400 µL
Hepes Buffer (25 mM), pH = 7.4	10 mL
Total Volume	10.53 mL

SAMPLE EXPERIMENTAL PROTOCOL

1. Mix equal volume of the alpha-thrombin standards or samples with 2X fluorescent alpha-thrombin substrate assay solution and incubate the solutions at room temperature for at least 1 hour.
2. Monitor the fluorescence using fluorescent microplate readers at 360/470 nm.

EXAMPLE DATA ANALYSIS AND FIGURES

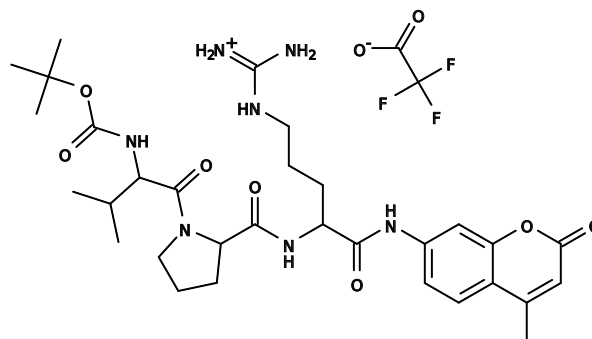


Figure 1. Chemical structure for BOC-Val-Pro-Arg-AMC

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.