

PRODUCT INFORMATION



Z-FR-AMC (trifluoroacetate salt)

Item No. 34045

Formal Name: N-[(phenylmethoxy)carbonyl]-L-phenylalanyl-N-(4-methyl-2-oxo-2H-1-benzopyran-7-yl)-L-argininamide, trifluoroacetate salt

Synonyms: Z-FR-7-amino-4-Methylcoumarin, Z-Phe-Arg-AMC

MF: $C_{33}H_{36}N_6O_6 \cdot XCF_3COOH$

FW: 612.7

Purity: $\geq 98\%$

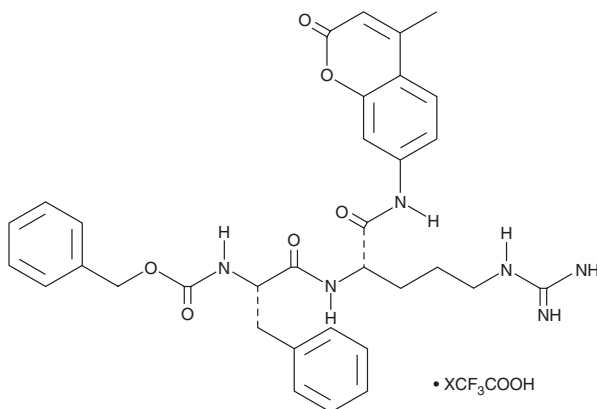
UV/Vis.: λ_{max} : 325 nm

Ex./Em. Max 340-360/440-460 nm

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Z-FR-AMC (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Z-FR-AMC (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Z-FR-AMC (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Z-FR-AMC (trifluoroacetate salt) in these solvents is approximately 1 mg/ml.

Description

Z-FR-AMC is a fluorogenic substrate for cathepsin L.¹ Upon enzymatic cleavage by cathepsin L, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify cathepsin L activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

Reference

1. Carmona, C., Dowd, A.J., Smith, A.M., *et al.* Cathepsin L proteinase secreted by *Fasciola hepatica* in vitro prevents antibody-mediated eosinophil attachment to newly excysted juveniles. *Mol. Biochem. Parasitol.* **62**(1), 9-17 (1993).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/13/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM