

PRODUCT INFORMATION



Z-LRGG-AMC (trifluoroacetate salt)

Item No. 26641

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

Synonyms: Z-Leu-Arg-Gly-Gly-AMC, Z-Leu-Arg-Gly-Gly-7-amino-4-methylcoumarin

MF: C₃₄H₄₄N₈O₈ • XCF₃COOH

FW: 692.8

Purity: ≥95%

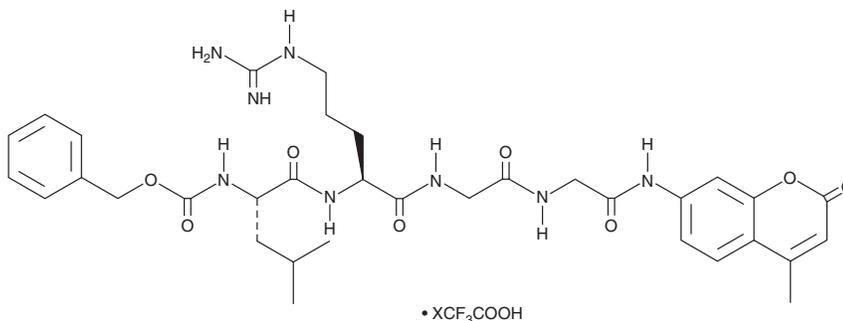
Ex./Em. Max: 340-360/440-460 nm

UV/Vis.: λ_{max}: 327 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Z-LRGG-AMC (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Z-LRGG-AMC (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Z-LRGG-AMC (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of Z-LRGG-AMC (trifluoroacetate salt) in these solvents is approximately 16 and 20 mg/ml, respectively.

Z-LRGG-AMC (trifluoroacetate salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Z-LRGG-AMC (trifluoroacetate salt) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Z-LRGG-AMC (trifluoroacetate salt) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

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

Reference

1. Stein, R.L., Chen, Z., and Melandri, F. Kinetic studies of isopeptidase T: Modulation of peptidase activity by ubiquitin. *Biochemistry* **34**(39), 12616-12623 (1995).

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, 11/04/2022

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