

PRODUCT INFORMATION



Ac-QQQ-AMC (hydrochloride)

Item No. 27711

Formal Name: (S)-2-acetamido-N¹-((S)-5-amino-1-(((S)-5-amino-1-((4-methyl-2-oxo-2H-chromen-7-yl)amino)-1,5-dioxopentan-2-yl)amino)-1,5-dioxopentan-2-yl)pentanediamide, hydrochloride

Synonyms: Ac-Gln-Gln-Gln-AMC,
Ac-Gln-Gln-Gln-7-amino-4-methylcoumarin,
Ac-L-Gln-L-Gln-L-Gln-AMC

MF: C₂₇H₃₅N₇O₉ • XHCl

FW: 601.6

Purity: ≥95%

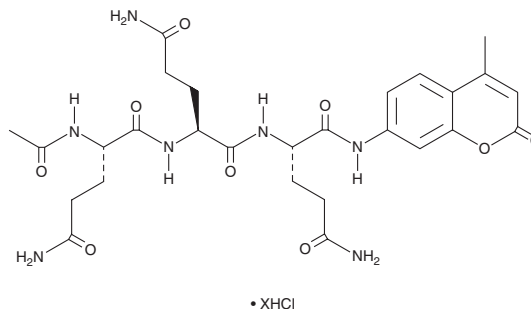
Ex./Em. Max: 340-360/440-460 nm

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

Ac-QQQ-AMC (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the Ac-QQQ-AMC (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Ac-QQQ-AMC (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Ac-QQQ-AMC (hydrochloride) in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

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

Description

Ac-QQQ-AMC is a fluorogenic probe. L-Glutamine tripeptide (QQQ) is a substrate for peptidoglutaminase I and II from the bacterium *B. circulans*.¹ AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

Reference

1. Kikuchi, M. A new enzymic method for determination of glutamine in proteins or peptides using peptidoglutaminases. *Anal. Biochem.* **59(1)**, 83-90 (1974).

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, 10/26/2022

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