

# PRODUCT INFORMATION



## Ac-RGK(Ac)-AMC (acetate)

Item No. 30798

**Formal Name:** (S)-6-acetamido-2-(2-((S)-2-acetamido-5-guanidinopentanamido)acetamido)-N-(4-methyl-2-oxo-2H-chromen-7-yl)hexanamide, acetate

**Synonym:** Ac-Arg-Gly-Lys(Ac)-AMC

**MF:** C<sub>28</sub>H<sub>40</sub>N<sub>8</sub>O<sub>7</sub> • XC<sub>2</sub>H<sub>4</sub>O<sub>2</sub>

**FW:** 600.7

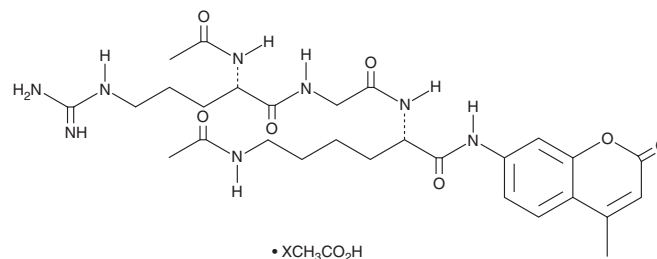
**Purity:** ≥98%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

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



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

### Laboratory Procedures

Ac-RGK(Ac)-AMC (acetate) is supplied as a solid. A stock solution may be made by dissolving the Ac-RGK(Ac)-AMC (acetate) in the solvent of choice, which should be purged with an inert gas. Ac-RGK(Ac)-AMC (acetate) is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of Ac-RGK(Ac)-AMC (acetate) can be prepared by directly dissolving the solid in aqueous buffers. Ac-RGK(Ac)-AMC (acetate) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Ac-RGK(Ac)-AMC is a fluorogenic substrate for histone deacetylases (HDACs).<sup>1</sup> Upon enzymatic cleavage by HDACs, 7-amino-4-methylcoumarin (AMC) is released and its fluorescence can be used to quantify HDAC activity. AMC displays excitation/emission maxima of 340-360/440-460 nm, respectively.

### Reference

1. Wegener, D., Wirsching, F., Riester, D., *et al.* A fluorogenic histone deacetylase assay well suited for high-throughput activity screening. *Chem. Biol.* **10**(1), 61-68 (2003).

#### 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, 05/22/2025

#### 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