

Ubiquitin-AMC

Ubiquitin substrate



Cat. No. 60-0116-050
Lot. No. 30399

Quantity: 50 µg
Storage: -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS

CERTIFICATE OF ANALYSIS Page 1 of 1

Background

In addition to fusion proteins, ubiquitin derivatives conjugated with a fluorophore have been reported as substrates for biochemical DUB assays. A frequently used coumarin-based substrate is ubiquitin-7-amido-4-methylcoumarin (Ub-AMC). DUBs catalyze the release of the AMC moiety, which is directly attached to the C-terminus of ubiquitin, and liberation of the fluorophore results in de-quenching of the fluorescent signal (Hassiepen *et al.*, 2007). The excitation/emission range of this fluorophore is 380nm/460nm respectively. The use of this substrate for determining steady-state kinetic parameters in a number of DUB assays was first described by Dang *et al.* (1998).

References:

Dang LC, Melandri FD and Stein RL (1998) Kinetic and mechanistic studies on the hydrolysis of ubiquitin C-terminal 7-amido-4-methylcoumarin by deubiquitinating enzymes. *Biochemistry* 37, 1868-1879.

Hassiepen U, Eidhoff U, Meder G, Bulber JF, Hein A, Bodendorf U, *et al.* (2007) A sensitive fluorescence intensity assay for deubiquitinating proteases using ubiquitin-rhodamine110-glycine as substrate. *Anal Biochem* 371, 201-207.

Physical Characteristics

Species: human

Source: synthetic

Quantity: 50 µg

Concentration: 2 mg/ml

Formulation: DMSO

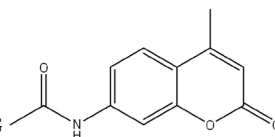
Molecular Weight: 8.72 kDa

Purity: >98% by InstantBlue™ SDS-PAGE

Stability/Storage: 12 months at -70°C; aliquot as required

Protein Sequence:

MQIFVKLTLTGKTTITLEVEPSDTIEN
VKAKIQDKEGIPPDQQLIFAGKQL
EDGRTLSDYNIQKESTLHLVLRLRGG

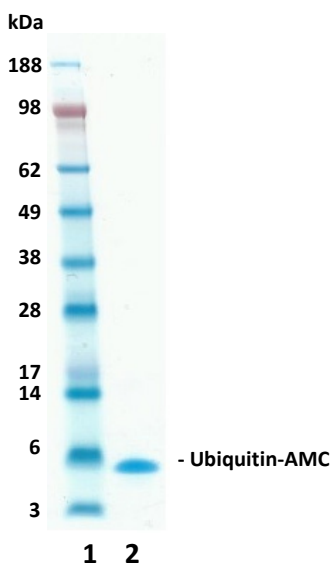


Ubiquitin (amino acid residues 1-76) C-terminally tagged with AMC (7-amino-4-methylcoumarin)
Accession number: P62987

Quality Assurance

Purity:

4-12% gradient SDS-PAGE
InstantBlue™ staining
Lane 1: MW markers
Lane 2: 1 µg Ubiquitin-AMC



Protein Identification:

Confirmed by mass spectrometry.

Activity Assay:

The activity of Ubiquitin-AMC (7-amido-4-methylcoumarin) was validated by determining the increase in fluorescence at 460nm (Excitation 380nm). Increased fluorescence is a result of the enzyme catalysed cleavage between the C-terminal Glycine and AMC, creating Ubiquitin and de-quenched AMC. UCHL3 (deubiquitylase) was incubated with Ubiquitin-AMC and the fluorescence was measured at four time points (0min, 30min, 60min and 90min).



www.ubiquigent.com
Dundee, Scotland, UK

ORDERS / SALES SUPPORT

International: +1-617-245-0003
US Toll-Free: 1-888-4E1E2E3 (1-888-431-3233)
Email: sales.support@ubiquigent.com

UK HQ and TECHNICAL SUPPORT

International: +44 (0) 1382 381147 (9AM-5PM UTC)
US/Canada: +1-617-245-0020 (9AM-5PM UTC)
Email: tech.support@ubiquigent.com

Email services@ubiquigent.com for enquiries regarding compound profiling and/or custom assay development services.

© Ubiquigent 2015. Unless otherwise noted, Ubiquigent, Ubiquigent logo and all other trademarks are the property of Ubiquigent, Ltd.

Limited Terms of Use: For research use only. Not for use in humans or for diagnostics. Not for distribution or resale in any form, modification or derivative OR for use in providing services to a third party (e.g. screening or profiling) without the written permission of Ubiquigent, Ltd.

Lot-specific COA version tracker: v1.0.0