Ubiquitin-Rhodamine 110

Ubiquitin substrate

Cat. No. 60-0117-050

Lot. No. 30371

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. Ubiquitin-Rhodamine 110 (Ub-Rho110-G) is a fluorogenic rhodamine-based substrate. While the disubstituted rhodamine moiety in Ub-Rho110-G is essentially non-fluorescent, cleavage results in a mono-substituted rhodamine, Rho110-G, which exhibits intense fluorescence when excited at 485 nm (Hassiepen et al., 2007). The rhodamine fluorophore exhibits optical properties more appropriate - than Ubiquitin-AMC - for compound screening and profiling. The risk of artifacts in screens due to autofluorescence of compounds is substantially reduced as the rhodamine 110 fluorophore has excitation and emission wavelengths of 485nm and 535nm respectively (Hassiepen et al., 2007).

References:

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

50 µg

-70°C

Species: human

Quantity:

Storage:

Source: synthetic

Quantity: 50 µg

Concentration: 2 mg/ml

Formulation: DMSO

Molecular Weight: 8.93 kDa

Purity: >95% by InstantBlue™ SDS-PAGE

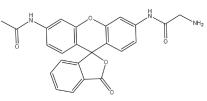
Stability/Storage: 12 months at -70°C;

aliquot as required

Protein Sequence:

MQIFVKTLTGKTITLEVEPSDTIEN VKAKIQDKEGIPPDQQRLIFAGKQL EDGRTLSDYNIOKESTLHLVLRLRGG

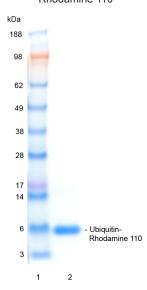
Ubiquitin (amino acid residues 1-76) C-terminally tagged with Rhodamine 110 Accession number: P62987



Quality Assurance

Purity:

4-12% gradient SDS-PAGE InstantBlue™ staining Lane 1: MW markers Lane 2: 1 µg Ubiquitin-Rhodamine 110



Protein Identification:

Confirmed by mass spectrometry.

Activity Assay:

The activity of Ubiquitin-Rhodamine 110 was validated by determining the increase in fluorescence at 535nm (Excitation 485nm) measured as a result of the enzyme catalysed cleavage at the amide bond between the C-terminal Glycine and Rhodamine, generating Ubiquitin and dequenched Rhodamine 110-Glycine. UCHL3 (deubiquitylase) was incubated with Ubiquitin-Rhodamine 110 and the fluorescence was measured at four time points (0min, 30min, 60min and 90min).



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

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

© **Ubiquigent 2014**. 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