

Ubiquitin-Rhodamine

High purity industry standard substrate for deubiquitinating enzymes (DUBs).

Ubiquitin-Rhodamine is prepared by the C-terminal derivatization of ubiquitin with modified rhodamine-110 and has been shown to be a useful and sensitive fluorogenic substrate for wide range of deubiquitinating enzymes (DUBs), including ubiquitin C-terminal hydrolases (UCHs) and ubiquitin specific proteases (USPs). Ubiquitin-rhodamine is particularly useful for studying deubiquitinating activity where detection sensitivity or continuous monitoring of activity is essential. The long excitation and emission wavelengths of rhodamine 110 (Excitation 485nm / Emission 520nm) avoid interference often caused at shorter wavelengths by small molecules and tissue culture components.

Ordering Information

[Order Online »](#)

BML-SE761-0025	25µg
----------------	------

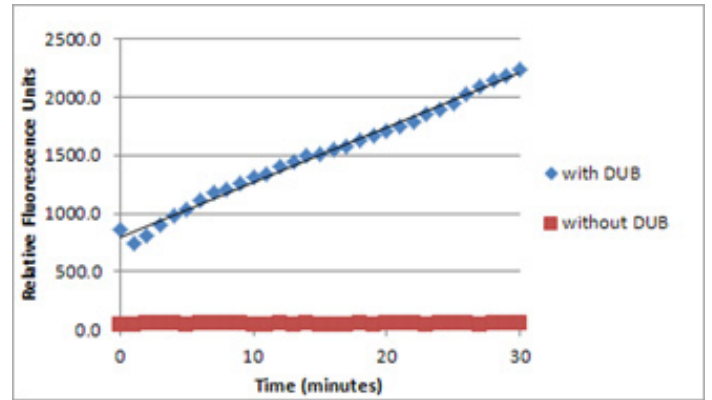
Manuals, SDS & CofA

[View Online »](#)



HPLC analysis of BML-SE761.

HPLC analysis of BML-SE761.



DUB Assay: 50nM USP2 (Prod. No. BML-UW9850); 500nM Ubiquitin-Rhodamine; 50mM HEPES pH7.8, 0.5mM EDTA, 1mM DTT, and ovalbumin 1mg/ml.

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least one year after receipt when stored at -80°C. After reconstitution, prepare aliquots and store at -80°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	-80°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application Notes	Can be used for the following applications: 1. Substrate for deubiquitylating enzyme activity assays. 2. Identification/confirmation of enzyme deubiquitylation activity. 3. Investigation of deconjugating enzyme substrate specificity in comparison with alternative UBL substrates.
Emission Maximum	Excitation 485nm / Emission 520nm
Formulation	Liquid. In dimethylsulfoxide.
MW	~8.9kDa
Purity	≥95% (HPLC)
Purity Detail	Please note: considerable efforts are taken to ensure that ubiquitin-rhodamine is homogeneous and not contaminated with ubiquitin and ubiquitin deletions, which are common in material obtained from other sources. The chromatogram below demonstrates the homogeneity of Prod. No. BML-SE761.
Sequence	Accession number: P62988; length: 76 amino acid residues.
Source	Produced in <i>E. coli</i> and covalently attached to modified rhodamine-110.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZOLIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
341 Benilouba
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 1 20 76 61 61
info-fr@enzolifesciences.com

Germany
Phone: +49 221 54 15 52
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

P0CG47 (UBB), P0CG48 (UBC), P62979 (RPS27A), P62987 (UBA52)