

Di-ubiquitin (K⁶³-linked)

Polyubiquitin chains, linked through specific lysine residues, are useful tools for investigating, amongst other things, the specificity and reactivity of deubiquitinating enzymes (DUBs) and the recognition and interaction of polyubiquitin modified proteins with by ubiquitin binding domain (UBDs) containing proteins.

Citations: 3

[View Online »](#)

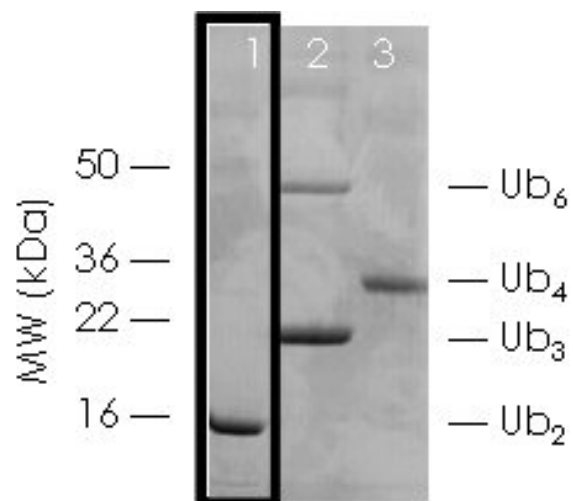
Ordering Information

[Order Online »](#)

BML-UW0730-0050	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE of di-ubiquitin Ub₂, K⁶³-linked (Lane 1: Prod. No. BML-UW0730, 5μg), in comparison with Ub₃, K⁶³-linked (Lane 2: Prod. No. [BML-UW0745](#), 5μg) and Ub₄, K⁶³-linked (Lane 3: Prod. No. BML-UW0715

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 12 months after receipt when stored at -20°C. It is recommended that the product be kept at a moderately high protein concentration (>1mg/mL) and in the presence of buffer. If a more dilute reagent is required, the inclusion of an inert carrier protein (e.g. ovalbumin at 0.5-1.0mg/mL) in the storage buffer is recommended in order to minimise loss of the reagent due to absorption.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ub2
Application Notes	<p>For use in deubiquitinating enzyme assays and polyubiquitin binding studies.</p> <p>Suggested uses:</p> <ol style="list-style-type: none">1. Deubiquitinating enzyme substrates (general/linkage specific).2. Investigation of polyubiquitin chain recognition by and interaction with ubiquitin binding proteins.
Formulation	Liquid. In 20mM TRIS-HCl, pH 7.5, containing 0.15M NaCl, 1mM EDTA.
MW	~17kDa
Purity	≥95% (SDS-PAGE)
Source	Protein components produced in <i>E. coli</i> . Synthesised enzymatically <i>in vitro</i> .

Technical Info / Product Notes

K⁶³-linked di-ubiquitin (Ub₂) is made by incubating E1 and Ubc13-Mms2 with equal concentrations of ubiquitin capped at the future proximal chain terminus (Asp⁷⁷-ubiquitin) and ubiquitin capped at the future distal chain terminus (K⁴⁸R/K⁶³R-ubiquitin). The distal terminus of the resulting dimer is de-blocked by treatment with yeast ubiquitin hydrolase-1 (YUH-1) followed by purification by anion exchange chromatography.

UniProt ID

P0CG47 (UBB), P0CG48 (UBC)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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