Di-ubiquitin (K⁶³-linked)

Polyubiquitin chains, linked through specific lysine residues, are useful tools for investigating, amongst other things, the specificity and reactivity of deubiquitinylating 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

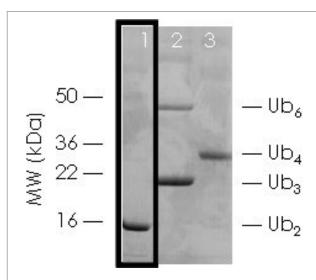
Order Online »

BML-UW0730-0050

50µg

Manuals, SDS & CofA

View Online »



SDS-PAGE of di-ubiquitin Ub $_2$, K 63 -linked (Lane 1: Prod. No. BML-UW0730, 5µg), in comparison with Ub $_3$, K 63 -linked (Lane 2: Prod. No. BML-UW0745, 5µg) and Ub $_4$, K 63 -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 For use in deubiquitinylating enzyme assays and

polyubiquitin binding studies.

Suggested uses:

Deubiquitinylating 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 *E. coli*. Synthesised

enzymatically in vitro.

Technical Info / Product Notes

 ${\rm K}^{63}\text{-linked di-ubiquitin (Ub}_2)$ is made by incubating E1 and Ubc13-Mms2 with equal concentrations of ubiquitin capped at the future proximal chain terminus (Asp⁷⁷ubiquitin) and ubiquitin capped at the future distal chain terminus ($K^{48}R/K^{63}R$ -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



info-