

Tetra-ubiquitin (K48-linked)

Ordering Information

[Order Online »](#)

ENZ-PRT266-0025	25µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Ub4

Formulation Liquid, containing 20mM Tris-HCl, pH 7.5, 0.15M NaCl, 1mM EDTA.

MW ~34kDa

Purity ≥90% (RP-HPLC)

Source Produced in *E. coli*.

Technical Info / Product Notes PLEASE NOTE: Tetra-ubiquitin may undergo complex formation to give higher molecular weight multimers upon heating and, consequently, the duration of any heating steps during use should be minimised. The exact nature of the multimeric complexes is not known at this time.

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)