

Background

Poly-ubiquitination is viewed as a posttranslational modification that controls protein stability or protein-protein interactions. Different poly-ubiquitination linkage determines the fate of modified protein(s). Increasing evidence suggests that polyubiquitin chains are generated by a ubiquitin E3 ligase complex, the linear ubiquitin chain assembly complex (LUBAC). The complex consists of 3 subunits, consisting of HOIL-1L, HOIP and SHARPIN. Linear polyubiquitin chains are cleaved by the deubiquitylases [USP2](#), CYLD, [USP5](#) and have been shown to bind to many ubiquitin binding domains (UBDs) including NEMO and Trabin-n (3xnfz). Recombinant tri-ubiquitin is expressed as a linear chain. amide linkages join the N- and C-terminus of each ubiquitin molecule to each other. This molecule is HIS-tagged at the N-terminus of the most distal ubiquitin.

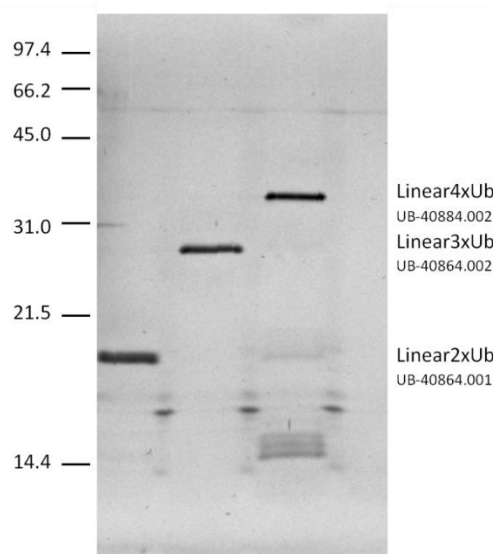
Application(s)

For use in deubiquitinating enzyme assays and polyubiquitin binding studies

Product Specifications

Affinity Tag	His6
Purity	≥ 90%
Molecular Weight	26.3 kDa
Quantity	100 µg
Species	Human
Expression System	<i>E. Coli</i>
Physical State	Liquid
Buffer	20mM Tris, pH 7.5, 0.15M NaCl, 1mM EDTA
Stability & Storage	1 year at -80° C. Avoid repeated freeze/thaw cycles

Product QC



Approximately 50ng of polyUb chains were subjected to SDS-PAGE analysis (15%) and visualized by silver staining.

References

1. Rittinger K, Ikeda F., et al., Open Biol. 2017;7(4):170026

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved

Linear tri-ubiquitin (Ub3)

Cat. # SI0103

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved