

# K27-linked di-ubiquitin (Ub2)

Cat. # SI2702

## Background

K27-linked diubiquitin (Ub2) is a diubiquitin molecule linked by a native isopeptide bond between the C-terminal glycine of the distal ubiquitin and the  $\epsilon$ -amino group of lysine 27 on the proximal ubiquitin. The product is generated using chemical ligation. Recent reports suggest that K27-linked chains play a role primarily in the intracellular innate immune response pathway, as well as in the DNA damage repair response. It is a useful substrate for identifying and characterizing deubiquitinating enzymes that specifically cleave the K27 linkage, and for structural and binding studies of ubiquitin chain recognition by ubiquitin-associated domains (UBA domains) or ubiquitin-interacting motifs (UIMs).

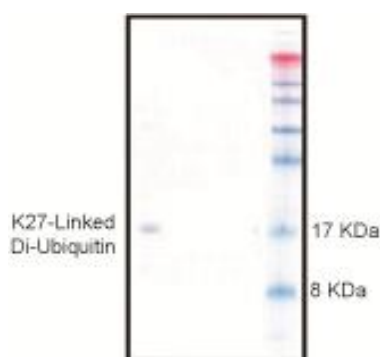
**Alternate Names** DiUbiquitin, Ub2

**Application(s)** Investigation of DUB linkage specificity.

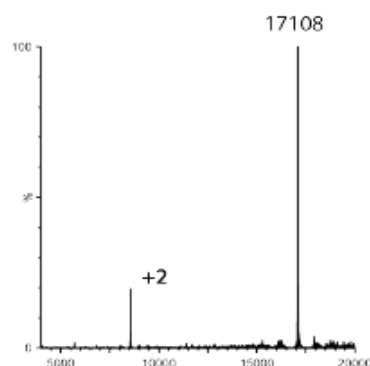
## Product Specifications

|                          |  |
|--------------------------|--|
| <b>Tag</b>               | None   |
| <b>Purity</b>            | $\geq 95\%$ by RP-HPLC                         |
| <b>Molecular Weight</b>  | 17108 Da by MS (calculated 17112 Da)           |
| <b>Quantity</b>          | 25 $\mu$ g                                     |
| <b>Species</b>           | Human  |
| <b>Expression System</b> | <i>E. Coli</i>                                 |
| <b>Physical State</b>    | Liquid at 0.5 mg/ml                            |
| <b>Buffer</b>            | 20 mM Tris-HCl, pH 7.5, 0.15 M NaCl, 1 mM EDTA |
| <b>Storage</b>           | -80°C. Avoid repeated freeze/thaw cycles       |

## Product QC



**SDS-PAGE Analysis**



**Mass Spectrum**

## References

1. van Tilburg GBA, et al., Cell Chem Biol. 2021;28(2):191-201.

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