

## K63-Linked Di-Ubiquitin (Phosphorylated)

Cat. # SI6302P

### Background:

Ubiquitin Chains are essential components in the study of protein degradation pathways, protein trafficking, and cellular signaling processes. These polymeric chains of ubiquitin molecules play critical roles in regulating protein stability, localization, and activity. K63 chains traditionally play a role in intracellular signalling, trafficking, and autophagy. Phospho-ubiquitin chains represent a specialized class of polyubiquitin characterized by phosphorylation at Serine 65 and play a central role in mitophagy signaling pathways.

K63 Di-Ubiquitin (phosphorylated) is a dimeric chain of wild-type ubiquitin, wherein ubiquitin monomers are linked together via an isopeptide bond between Lysine 63 and the C-terminal Glycine. The chains are then enzymatically phosphorylated at the Ser65 position.

### Application:

- Investigation of phosphoubiquitin chain specificity and selectivity
- Studies on the role of phosphoubiquitin chains in protein degradation pathways (e.g., proteasomal and autophagic degradation)
- Analysis of phosphoubiquitin-mediated signaling pathways and cellular responses
- Structural studies to elucidate the architecture and dynamics of phosphoubiquitin chains
- Screening assays to identify modulators of phosphoubiquitin chain assembly and disassembly processes

### Product Information

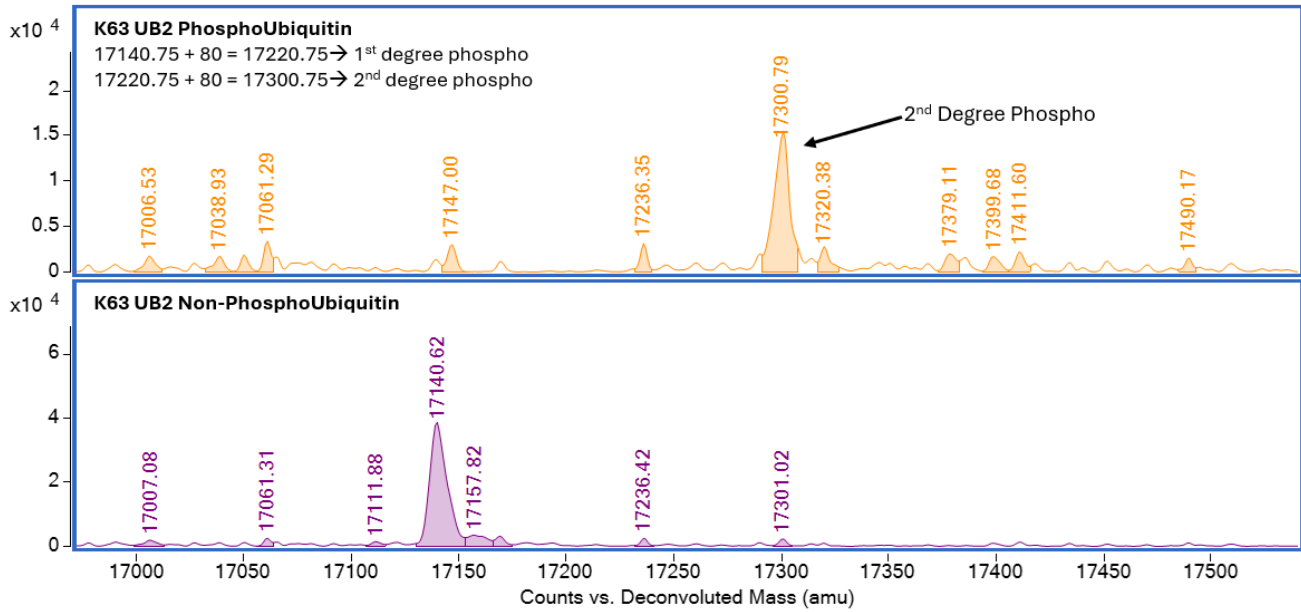
<b>Purity:</b>	≥ 95% by HPLC-MS
<b>Molecular Weight:</b>	17219-17299 Da
<b>Physical State:</b>	Liquid, 50 mM Tris, pH 7.5, 0.15 M NaCl
<b>Quantity:</b>	100 µg
<b>Solubility:</b>	>1 mg/mL
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles

### References

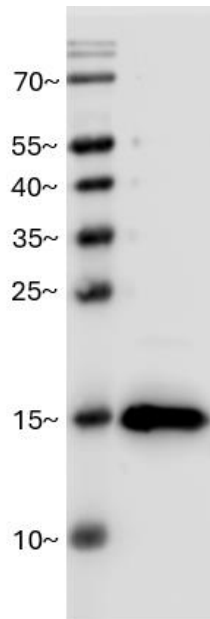
1. Madiraju, C.; Novack, J.P.; Red, J.C.; Matsuzawa, S-I. K63 Ubiquitination in Immune Signaling. *Trends Immunol.* **2022**, *43*, 148-162.
2. Cao, L.; Liu, X.; Zheng, B.; Xing, C.; Liu, J. Role of K63-linked Ubiquitination in Cancer. *Cell Death Discov.* **2022**, *8*, 410.
3. Swatek, K.N. & Komander, D. Ubiquitin Modifications. *Cell Res.* **2016**, *26*, 399-422.
4. Yau, R. & Rape, M. The increasing complexity of the ubiquitin code. *Nature Cell. Bio.* **2016**, *18*, 579-586.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2009 LifeSensors, Inc. All Rights Reserved

Data



Deconvoluted Mass Spectrum



Western Blot Analysis (100 ng)

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
 Copyright © 2009 LifeSensors, Inc. All Rights Reserved