

DUB Cocktail

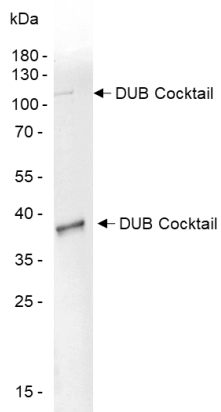
Cat. # DB599

Background This DUB cocktail contains an optimized combination of DUBs to be used as part of the [UbiTest](#) assay. Prior experiments with single broad-spectrum DUBs did not completely cleave all ubiquitin chains. This cocktail was designed for complete DUB digestion as part of any assay.

Product Information

Purity	≥ 95% by SDS-PAGE
Molecular Weight	Varies
Quantity	25 µg
Physical State	Liquid
Source	Human Recombinant
Tag	His6-SUMO
Activity	This enzyme is active in the UbiTest assay.
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified DUB Cocktail. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.

References

1. Lange, S., et al., Mol Cell., 2022. 82(1):15-29.
2. Mevissen, T., and Komander, D., Annu Rev Biochem., 2017. 86:159-192.

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