

OTUD7B (Cezanne)

Cat. # DB208

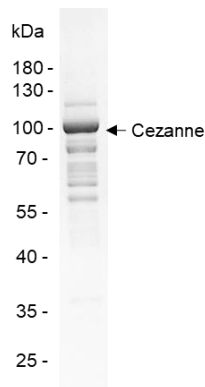
Background Cezanne (OTUD7B) is a deubiquitinase that preferentially cleaves K11 ubiquitin linkages. This enzyme was also shown to cleave K48- and K63-linked ubiquitin chains *in vitro*. It is known to be a negative regulator of NF- κ B signaling.

Alternate names OTU domain-containing protein 7B, ZA20D1

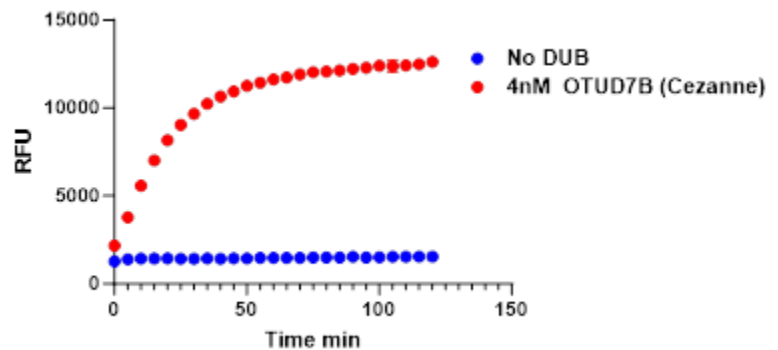
Product Information

Molecular Weight	103 kDa
Quantity	25 μ g
Purity	\geq 80% by SDS-PAGE
Physical State	Liquid
Source	Human Recombinant
Tag	His6-SUMO
Storage	-80 $^{\circ}$ C. Avoid repeated freeze/thaw cycles

Product QC



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



Activity Assay of Cezanne. 4 nM Cezanne was tested in Ub-Rh110 assay showing robust DUB activity.

References

1. Chen, X., et al., Am J Cancer Res., 2020. 10(12):4342-4356.
2. Wang, Y., et al., Cell Death Discov., 2023. 9(1):302.

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