

OTUD6B

Cat. # DB205

Background: OTUD6B is a deubiquitylase belonging to the OTU (ovarian tumor domain) family of DUBs. OTUD6B regulates the proliferation of B lymphocytes, in part by controlling the cell cycle through proteins such as cyclin D2 [1]. The ability of OTUD6B to regulate cell proliferation has been implicated in connection with non-small cell lung cancer (NSCLC) [2].

Alternate names: Human deubiquitinating enzyme ovarian tumor domain-containing 6B, DUBA5, OTUD-6B, OTUD6B (GenBank Accession No. NM_016023).

Product Information

Accession No.	NM_016023
Molecular Weight:	39 kDa
Quantity:	50 µg
Physical State:	Liquid, 25 mM Tris-HCl, pH 7.5, 150 mM NaCl, 5 mM DTT, 10% glycerol
Source:	Human Recombinant
Tag:	N-terminal His6-Flag tag
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

1. Xu, Y., Zheng, Y., Zhu, Y., Kong, X., and Hu, L., Evidence for OTUD-6B Participation in B Lymphocytes Cell Cycle After Cytokine Stimulation. *PLoS One*. **6**: e14514 (2011).
2. Sobol, A., Askonas, C., Alani, S., Weber, M.J., Anathanarayanan, V., Osipo, C., and Bocchetta, M., Deubiquitinase OTUD6B Isoforms are Important Regulators of Growth and Proliferation. *Mol Cancer Res*. **15**:117-127 (2017).

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