

USP14 (Ubiquitin Specific Protease 14)

Cat. # DB505

-
- Background:** USP14 is an ubiquitin specific protease that associates with the proteasome. This association has been shown to enhance the enzymatic activity of the proteasome^{3,4}. Recent studies have demonstrated that USP14 expression in certain cancers correlates with metastasis². Using a chemical probe researchers have shown that USP14 can interact with ISG15, a ubiquitin-like protein¹.
- Alternate names:** Deubiquitinating enzyme 14, TGT, Ubiquitin carboxyl-terminal hydrolase 14, Ubiquitin-specific-processing protease 14, Ubiquitin thioesterase 14
-

Product Information

Accession No.:	P54578
Molecular Weight:	68kDa
Quantity:	25µg
Physical State:	Liquid
Source:	Human, recombinant
Tag:	His ₆ SUMO
Activity:	This enzyme cleaves K48-linked ubiquitin chains in vitro.
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

1. Catic, A. et al. Screen for ISG15-crossreactive deubiquitinases. *PLoS ONE* **2**, e679 (2007).
2. Shinji, S. et al. Ubiquitin-specific protease 14 expression in colorectal cancer is associated with liver and lymph node metastases. *Oncol Rep* **15**, 539-43 (2006).
3. Hu, M. et al. Structure and mechanisms of the proteasome-associated deubiquitinating enzyme USP14. *EMBO J* **24**, 3747-3756 (2005).
4. Borodovsky, A. et al. A novel active site-directed probe specific for deubiquitylating enzymes reveals proteasome association of USP14. *EMBO J* **20**, 5187-5196 (2001).

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