

# USP2 (catalytic domain) (rat), (recombinant) (untagged)

Ubiquitin-specific proteases (USPs) represent the most widespread and represented deubiquitylating enzymes across evolution. USPs tend to release ubiquitin from a conjugated protein. They display similar catalytic domains but divergent N-terminal and occasionally C-terminal extensions, which are thought to function in substrate recognition, subcellular localization, and protein-protein interactions.

USP2 is a ubiquitin-specific protease thought to function at the pre-proteasomal level by preventing the degradation of fatty acid synthase (FAS). FAS overexpression occurs in the majority of epithelial tumors, including prostate cancer, and protects cancer cells from apoptosis. Because USP2 is overexpressed in prostate cancer and its inhibition results in apoptosis of transformed prostatic cells, this isopeptidase is a potential target for induction of apoptosis in prostate cancer cells.

The full length USP2 protein is comprised of 618 amino acids with a predicted molecular mass of 69.4 kDa and contains the highly conserved Cys and His boxes present in all members of the UBP family of deubiquitylating enzymes. The N-terminal region (residues 271-618) of USP2 contains all the necessary residues for catalysis and exhibits deubiquitylating activity vs. ubiquitin-AMC (Prod. No. BML-SE211).

Citations: 6

[View Online »](#)

## Ordering Information

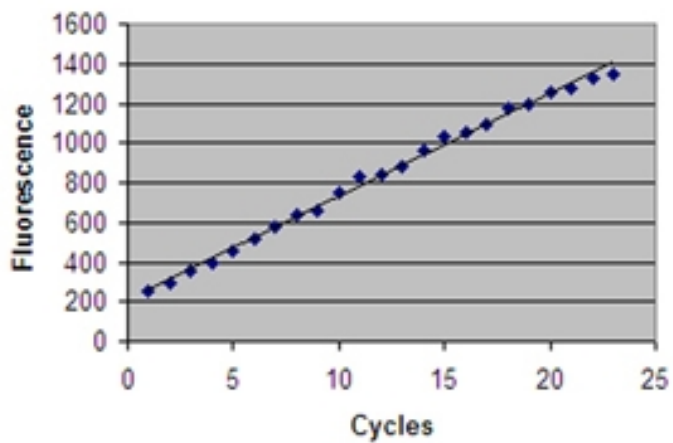
[Order Online »](#)

BML-UW9850-0100

100µg

## Manuals, SDS & CofA

[View Online »](#)



Example of assay result. DUB assay: 50 nM USP2 catalytic domain (BML-UW9850); 500nM Ubiquitin-AMC; 50mM HEPES, pH 7.8, 0.5mM EDTA, 1mM DTT



1 2

Coomassie stained gel of USP2 catalytic domain (UW9850). Lane 1: MW markers (top to bottom) 205, 116, 97, 84, 66, 55,45 and 36. Lane 2: UW9850 (2ug).

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

**Alternative Name** Ubiquitin-specific protease 2

### Application Notes

#### Uses:

Deubiquitinylation assays using small molecule substrates (e.g. Ub-AMC) or ubiquitin-protein conjugates.  
Studies of the interaction of USP2 with FAS1 and their role in prostate cancer.  
Development of inhibitors for USP2 deubiquitinating activity.

### Formulation

Liquid. In 10mM MES, 0.5mM DTT, 50mM TRIS-HCl, pH 7.5, containing 20% glycerol.

### MW

~40kDa

### Purity

≥95% (SDS-PAGE, following ion exchange chromatography)

### Source

Produced in *E. coli* BL21 (λDE3). Untagged USP2 catalytic domain.

### UniProt ID

Q5U349

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)