

Ubiquitin vinyl sulfone, (HA-tag)

For use in the detection and identification of deubiquitylating enzymes. HA-Ub-VS is a DUB active site directed probe generated by a chemical ligation method, that acts as a potent and irreversible inhibitor of DUBs through covalent modification of the active site and as a specific probe for enzymes with DUB activity. The HA peptide sequence (YPYDVPDYA), derived from the influenza hemagglutinin protein, facilitates sensitive identification or purification of HA-Ub-VS modified DUBs through recognition by anti-HA antibodies and/or anti-HA-agarose.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-UW0155-0025	25µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 6 months after receipt when stored at -80°C. Once reconstituted store HA-Ub-VS solution at -80°C, avoid multiple freeze/thaw cycles and use within 1 month.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	-80°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	HA-Ub-VS
Application Notes	<p>Suggested uses:</p> <ol style="list-style-type: none">1. Deubiquitinating enzyme active site-directed probes for detection of novel DUBs.2. Determination of deubiquitinating enzyme activity of putative DUBs.3. Identification of DUBs from biological samples via immunoprecipitation and proteomic methods.4. Characterization of deconjugating enzymes for Ub/Ubl specificity.
Formulation	Lyophilized.
MW	~9.8kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with aqueous buffer up to 5mg/ml (~500µM).
Source	Synthetic.

Technical Info / Product Notes

Typical assay set-up: HA-Ub-VS probe concentration: 0.5-5.0µM. Purified enzyme concentration: 0.1-1.0µM; Buffer: 50mM TRIS, pH7.5, 150mM NaCl, 1mM DTT. Incubate at 37°C for 30 minutes. Modification of DUB enzymes by site-directed HA-Ub-VS probes detected by Western blotting with either HA antibody (general reactivity) or an antibody specific to particular DUB of interest. **Note:** When probing complex solutions such as cell lysates the optimal amount of total sample protein required will depend on the specific lysate to be analyzed and must therefore be determined for the individual lysate of interest.

UniProt ID

P0CG47 (UBB), P0CG48 (UBC), P62979 (RPS27A), P62987 (UBA52)

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-
be@enzolifesciences.com](mailto:info-be@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)