

# SUMO-QAPTURE-T®

## kit

The only commercially available kit for isolation and enrichment of SUMOylated proteins.

The SUMO-QAPTURE-T® kit is an efficient tool for the selective isolation of SUMOylated proteins. The kit facilitates the affinity purification of SUMOylated proteins from cell extracts and tissue lysates using a high-binding SIM-containing affinity matrix. Captured proteins are eluted under denaturing conditions followed by analysis by Western blotting, using the SUMO antibodies provided or antibodies to specific proteins of interest, or potential substrate identification by proteomic methods. SUMO-conjugate-containing samples are prepared in native form in the presence of protease inhibitors to prevent loss through the action of deSUMOylating enzymes. Optimization of binding permits complete isolation of the full range of SUMO-protein conjugates from a specific lysate.

Citations: 9

[View Online »](#)

### Ordering Information

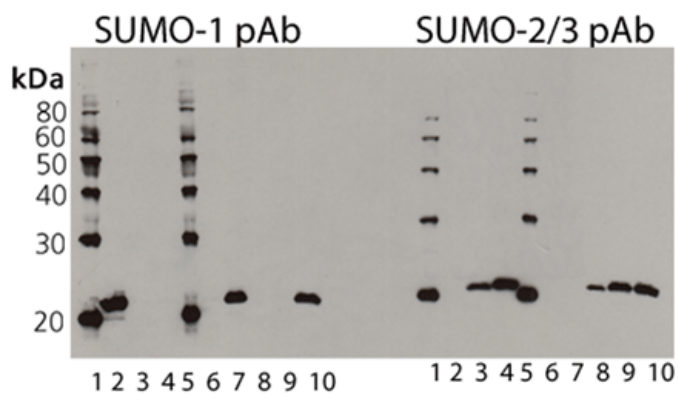
[Order Online »](#)

BML-UW1000A-0001	10 tests
------------------	----------

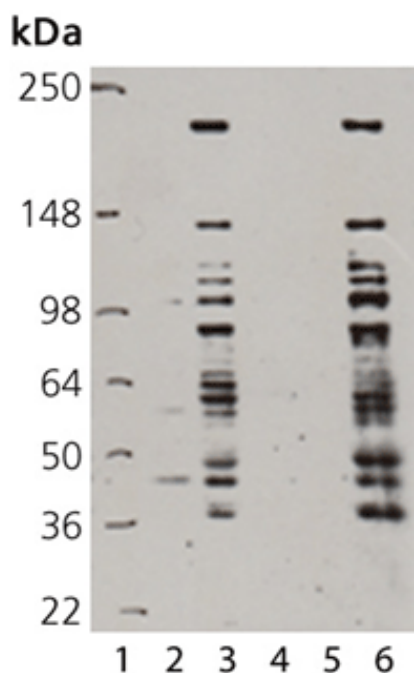
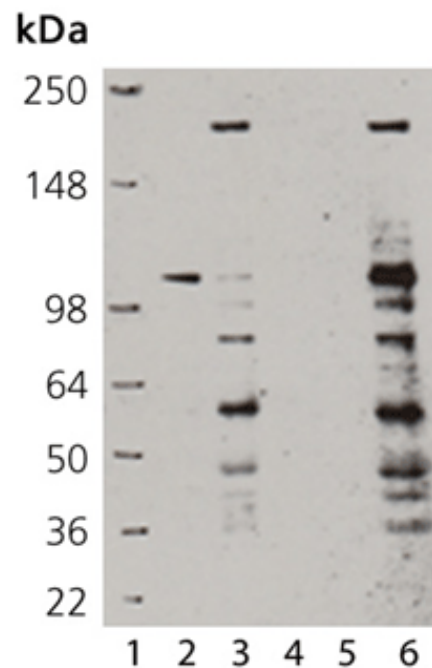
### Manuals, SDS & CofA

[View Online »](#)

- Fast, simple, and reliable assay
- High specificity, high through-put capacity
- Amenable to analysis via Western blotting or proteomic methods



Western blot analysis of SUMO-QAPTURE-T<sup>®</sup> binding of recombinant SUMO protein controls. 200 ng of each recombinant control were bound alone or simultaneously using the described protocol. Lane 1: MWM, Lane 2: 40 ng SUMO-1, Lane 3: 40 ng SUMO-2, Lane 4: 40 ng SUMO-3, Lane 5: MWM, Lane 6: SUMO 1/2/3 final wash fraction, Lane 7: SUMO-1 elution, Lane 8: SUMO-2 elution, Lane 9: SUMO-3 elution, Lane 10: SUMO 1/2/3 elution. Left probed with SUMO-1 pAb, Right probed with SUMO 2/3 pAb.



## Handling & Storage

Use/Stability	Unopened kit should be stored at -80°C upon receipt. After thawing, SUMO-QAPTURE-T® matrix should be stored at +4°C. Avoid a second freeze/thaw cycle. SUMO-1, SUMO-2 and SUMO-3 controls should be stored at -80°C. SUMO antibody solutions can be stored at -20°C. Avoid multiple freeze/thaw cycles of components to ensure stability and activity.
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

Application Notes	<p>For the isolation and enrichment of SUMOylated proteins.</p> <p>Suggested Applications:</p> <ol style="list-style-type: none"><li>1. Capture and isolation of SUMO-protein conjugates from specific cell/tissue lysates of interest with subsequent detection and analysis by Western blotting.</li><li>2. Identification of SUMO-modified protein substrates by proteomic analysis methods following release of free SUMOylated proteins in denatured form.</li><li>3. Selective purification/pull down of SUMOylated proteins from <i>in vitro</i> SUMOylation assays.</li></ol>
Contents	<p>SUMO-QAPTURE® matrix (BML-UW0980-0200)</p> <p>SUMO-1 Control (BML-UW9195-0125)</p> <p>SUMO-2 Control (BML-UW9205-0125)</p> <p>SUMO-3 Control (BML-UW9215-0125)</p> <p>SUMO antibody solutions includes,</p> <p>SUMO-1 polyclonal antibody (BML-PW0505A-0010)</p> <p>SUMO-2 polyclonal antibody (BML-PW0510A-0010)</p>
Quantity	Sufficient for 10 binding assays.
Technical Info / Product Notes	The former SUMOylated-protein lysate control (Prod. No. BML-UW0130A) is available upon request.



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)