

# SUMOylation kit

Market leading kit for generation of SUMOylated proteins *in vitro*.

This kit provides a means of generating SUMOylated proteins *in vitro*, by covalent linkage of the carboxy-terminal of SUMO-1, -2 or -3 to specific lysine residues on the target protein *via* isopeptide bonds, using the SUMOylation enzyme cascade. A short sequence containing the consensus  $\Psi$ -K-X-D/E (where lysine is the amino acid modified,  $\Psi$  is a large hydrophobic residue and X is any amino acid residue) is thought to be necessary for this *in vitro* protein SUMOylation to occur, however SUMOylation has also been observed in cases where the consensus site is not conserved. A control target protein is provided together with all other necessary components. SUMO specific antibodies are provided for detection of SUMOylated proteins *via* SDS-PAGE and Western blotting. Provides sufficient material for 20 x 20 $\mu$ L reactions. Suggested uses: For SUMO-modification of specific proteins *in vitro*, To demonstrate that novel proteins are potential targets for SUMOylation under *in vitro* conditions, To generate substrates for deSUMOylating enzymes, such as SENP1 and SENP2, To test proteins for SUMO E3 ligase activity.

Citations: 59

[View Online »](#)

## Ordering Information

[Order Online »](#)

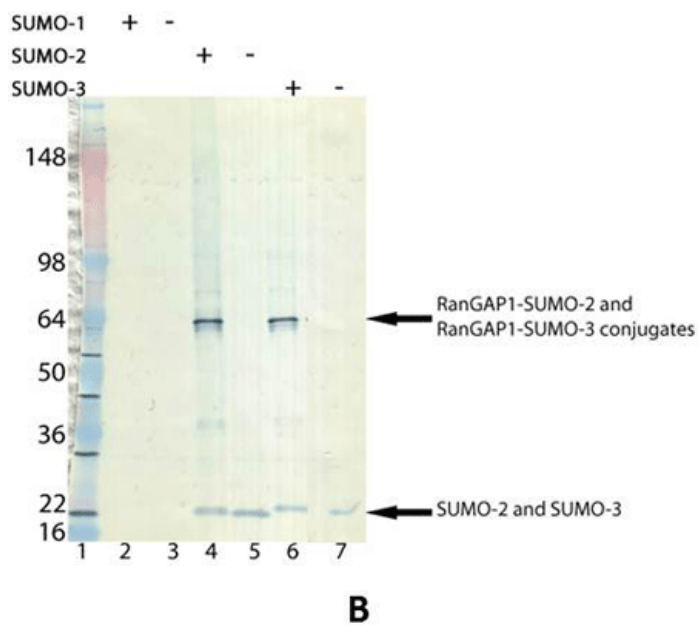
BML-UW8955-0001

20 tests

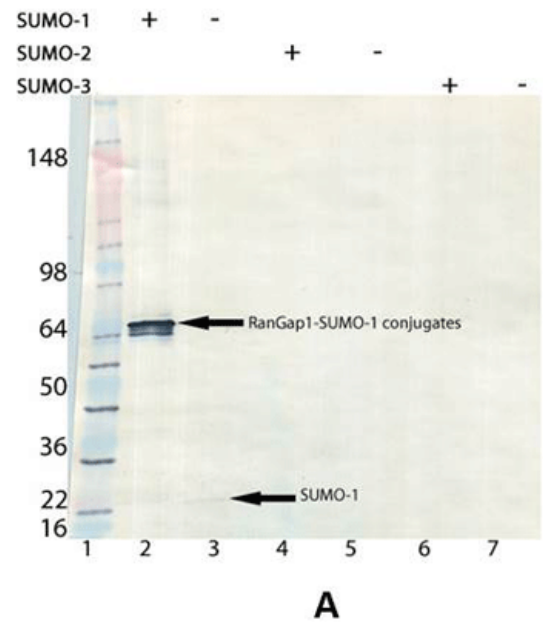
## Manuals, SDS & CofA

[View Online »](#)

- Fastest assay time on the market, just under 1.5 hours
- Simple 4 step assay: Mix, Incubate, Quench and Assay
- High specificity, high throughput capacity
- Amenable to analysis via Western blotting or proteomic methods



Western Blot of SUMOylation Assays for control RanGAP1 target protein. Assays set-up and run as described in "Assay Protocol". SUMOylated proteins were detected by Western Blotting as described in "Analysis by Western Blotting". B: detected by SUMO-2/3, pAb (BML-PW9465).



Western Blot of SUMOylation Assays for control RanGAP1 target protein. Assays set-up and run as described in "Assay Protocol". SUMOylated proteins were detected by Western Blotting as described in "Analysis by Western Blotting". A: detected by SUMO-1, pAb (BML-PW8330)

# Handling & Storage

Use/Stability	Unopened kit should be stored at -80°C to ensure stability and activity. After opening, The SUMO antibody solutions (BML-PW8330 and BML-PW9465) can be stored at -20°C or -80°C. Other components should be stored at -80°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

### Uses:

1. SUMO-modification of specific proteins *in vitro*. Allow investigation of the effect SUMOylation has on enzyme function, stabilisation, protein:protein interactions and, hence, its role in regulation of cellular processes, such as the p53 tumour repressor and NF- $\kappa$ B pathways.
2. Demonstrate novel proteins are potential targets for SUMOylation under *in vitro* conditions. Starting point for examining the role SUMOylation of a protein might play *in vivo*.
3. Generate substrates for deSUMOylating enzymes, such as SENP1 (Prod. No. BML-UW9760) and SENP2 (Prod. No. BML-UW9765).
4. Test proteins for SUMO E3 ligase activity: does it facilitate or enhance SUMOylation of specific target proteins, particularly under conditions/enzyme concentrations that more closely represent those *in vivo*.
5. Addition of known SUMO E3 ligase to facilitate/enhance target protein SUMOylation, particularly under conditions/enzyme concentrations that more closely represent those *in vivo* (e.g. RANBP2 [Prod. No. BML-UW9455], shown to be a ligase for SP100 SUMOylation).
6. SUMOylation of proteins in cell lysates or crude fractions/preparations to facilitate investigation of their role/function in complex solutions.
7. Demonstrate SUMOylation of known proteins in specific lysates (confirm with target protein specific antibodies).
8. Use of cell lysate or crude fractions/preparations as source of SUMO E3 ligases to facilitate SUMOylation of purified target proteins in the presence of SUMOylation kit components.

**Note:** Protocol provided for application 1. Assay set-up can be readily modified for alternative applications by inclusion, omission or substitution of specific components.

## Contents

### **20x SUMO Activating Enzyme Solution (SUMO E1):**

SUMO activating enzyme E1 (human), (recombinant), (Prod. No. BML-UW9330), 20 µl.

### **20x SUMO Conjugating Enzyme Solution (SUMO E2):**

Ubc9 (human), (recombinant) (untagged), (Prod. No. BML-UW9320), 20 µl.

### **20x SUMO Enzyme Solutions (SUMO-1, SUMO-2,**

**SUMO-3):** SUMO-1 (human), (recombinant) (His-tag), (Prod. No. BML-UW9195), 20 µl;

SUMO-2 (human), (recombinant) (His-tag), (Prod. No. BML-UW9205), 20 µl;

SUMO-3 (human), (recombinant) (His-tag), (Prod. No. BML-UW9215), 20 µl.

10x SUMOylation Buffer (Prod. No. BML-KW9890), 40 µl

### **20x Control RanGAP1 SUMOylation Target Protein**

**Solution:** RanGAP1 fragment (human), (recombinant) (GST-tag), (Prod. No. BML-UW9755), 20 µl.

**20x Mg-ATP Solution** (Prod. No. BML-EW9805), 25 µl

### **SUMO Antibody Solutions:**

SUMO-1 (human) (NT), polyclonal antibody, (Prod. No. BML-PW8330), 25 µl;

SUMO-2/3 (human) (NT), polyclonal antibody, (Prod. No. BML-PW9465), 25 µl.

## Quantity

Sufficient for 20 assays.

Last modified: September 3, 2025



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)