

# SUMO activating enzyme E1 (human), (recombinant)

SUMO conjugation to target proteins is mediated by a different, but analogous, pathway to ubiquitinylation. SUMO activating enzyme is a heterodimeric complex consisting of Aos1 and Uba2. Both subunits are well conserved from yeast to humans, with Aos2/SAE I being similar to the N-terminal half of the E1 enzyme for ubiquitin while Uba2/SAE II has similarity to the C-terminal half, and contains the active site cysteine residue required for formation of thioester bonds. However, Uba2 alone is not sufficient to catalyze SUMOylation. SUMO1 activating enzyme supports *in vitro* SUMOylation.

Citations: 5

[View Online »](#)

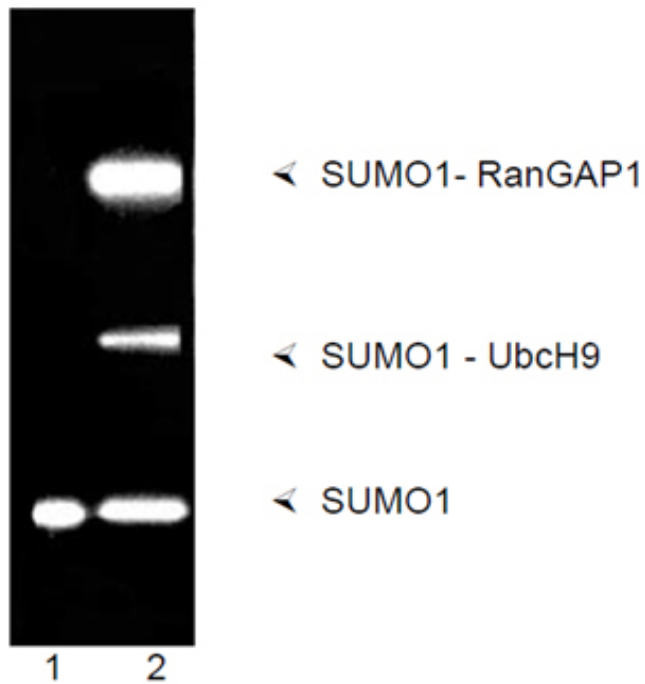
## Ordering Information

[Order Online »](#)

BML-UW9330-0025	25µg
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)



SUMOylation assay utilising SUMO-1, SUMO E1 (Prod. No. BML-UW9330), SUMO E2 (Prod. No. BML-UW9320) and RanGAP1 as target protein in presence (lane 2) and absence (lanes (1) of ATP after SDS-PAGE and blotting to PVDF with subsequent probing with Prod. No. BML-PW9460 (Anti-SUMO1).



SUMO E1 heterodimer by SDS-PAGE. Lane M: Markers, in kDa (from top to bottom), 205, 116, 97, 84, 66, 55, 45 and 36; Lane 1: SUMO E1 heterodimer, final product solution.

## Handling & Storage

Long Term Storage      -80°C

Shipping      Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Formulation**      Liquid. In 20mM HEPES, pH 8.0, containing 110mM potassium acetate, 2mM magnesium acetate, 0.5mM EGTA and 1mM DTT.

**MW**      ~38kDa and 90kDa

**Source**      Produced in *E. coli*. Full length heterodimeric human SUMO-activating enzyme.

**UniProt ID**      Q9UBE0 (subunit 1), O95605 (subunit 2)



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)