

# SUMO-4 (human), (recombinant) (His- tag)

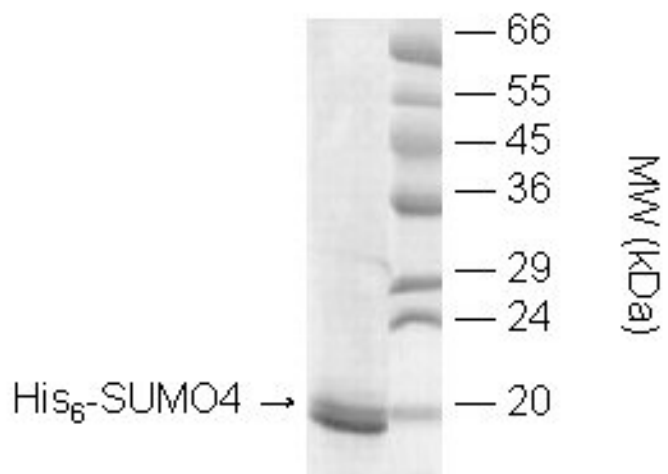
SUMO-4 is a member of the SUMO (small ubiquitin-related modifier) family of ubiquitin-like proteins involved in the regulation of the cellular functions of a wide variety of proteins, although there remains considerable debate as to whether this is a real and expressed gene product.

It has been reported that upon oxidative stress, SUMO-4 conjugates to various anti-oxidant enzymes, chaperones, and stress defence proteins, whereas under standard physiological conditions it is rapidly degraded. Subsequent proteomic analysis of SUMO-4 conjugates identified substrates implicated in the regulation of DNA repair and synthesis, RNA processing, protein degradation and glucose metabolism, suggesting a role for SUMO-4 in the regulation of intracellular stress.

The M55V substitution in SUMO-4 has been associated with type I diabetes.

**Manuals, SDS & CofA**

[View Online »](#)



SDS-PAGE image of His<sub>6</sub>-SUMO4 (1.5μg, BML-UW0905), coomassie stained.

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 12 months after receipt when stored at -80°C.
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
<b>Short Term Storage</b>	-80°C
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Small ubiquitin-related modifier 4
<b>Application Notes</b>	For use in SUMO-4 conjugation studies.
<b>Formulation</b>	Liquid. In 20mM HEPES, pH 8.0, containing 1mM DTT.
<b>MW</b>	~12.5kDa
<b>Purity</b>	≥90% (SDS-PAGE)
<b>Purity Detail</b>	Purified by multistep chromatography.
<b>Source</b>	Produced in <i>E. coli</i> .
<b>UniProt ID</b>	Q6EEV6

**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)