

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.

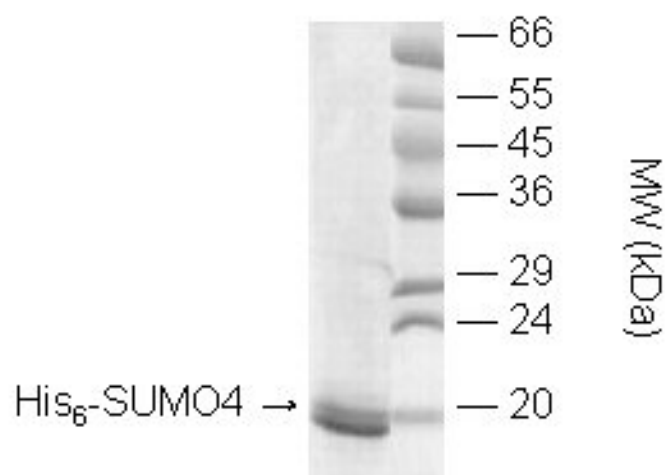
Ordering Information

[Order Online »](#)

BML-UW0905-0100	100µg
------------------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE image of His₆-SUMO4 (1.5μg, BML-UW0905), coomassie stained.

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

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