

Urm1 (human), (recombinant) (His- tag)

Urm1 is a member of the ubiquitin-like family of proteins and acts as a post-translational protein modifier in *Saccharomyces cerevisiae*. Simultaneous loss of Urm1p and Cla4p, a p21-activated kinase that functions in budding, is lethal, suggesting a role for the urmylation pathway in budding whilst additional results suggest an involvement in nutrient sensing. The first *in vivo* target for the urmylation pathway has been identified as the antioxidant protein Ahp1p. It has been suggested that the conjugation of Urm1p to Ahp1p could regulate the function of Ahp1p in antioxidant stress response in *Saccharomyces cerevisiae*.

Ordering Information

[Order Online »](#)

BML-UW9530-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

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, 50mM NaCl, 1mM DTT.
MW	16.3 kDa
Purity	≥95% (SDS-PAGE)
Source	Produced in <i>E. coli</i> .
UniProt ID	Q9BTM9

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)