

MetAP2 (human), (recombinant) (HA- tag)

MetAP2 is a metalloenzyme that catalyzes removal of the N-terminal methionine from proteins, with the physiologically relevant metal thought to be either cobalt or manganese. Since it was identified as the target of the anti-angiogenic natural product fumagilin (Prod. No. BML-CT100), it has been studied for its involvement in cancer, angiogenesis, inflammation, and rheumatoid arthritis. Because of differing substrate specificities and kinetics it is not desirable to substitute bacterial methionine aminopeptidase for mammalian MetAP2.

Ordering Information

[Order Online »](#)

BML-SE569-0010	10µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -80°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Methionine aminopeptidase 2, p67eIF2, Eukaryotic initiation factor 2-associated protein
Formulation	Liquid. In 20mM HEPES, pH 7.3, containing 0.1M KCl, 1mM CoCl ₂ , 1.5mM MgCl ₂ , 0.02% sodium azide and 10% glycerol.
MW	~70kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by chromatography
Source	Produced in insect cells. Recombinant, full-length (aa 1-478) methionine aminopeptidase 2 with an N-terminal hemagglutinin (HA) tag (YPYDVDPDYA), cloned from human cDNA. Produced in a baculovirus expression system.
UniProt ID	P50579



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)