

Cyclophilin A (human), (recombinant)

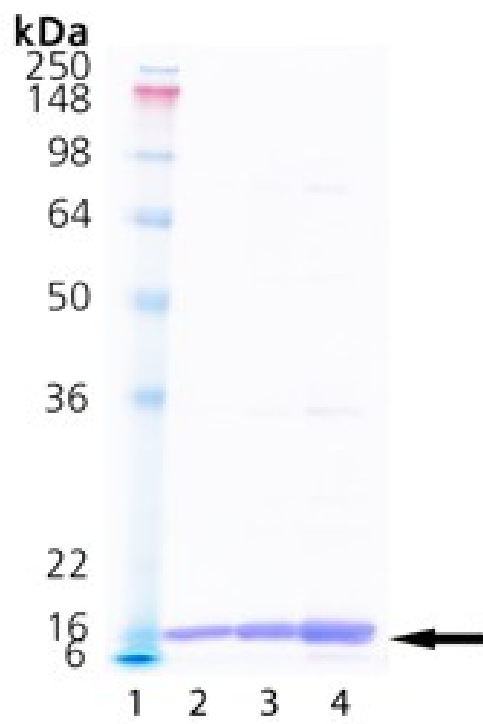
Ordering Information

[Order Online »](#)

BML-SE105-0100	100µg
-----------------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis. Lane 1: MW Marker. Lane 2: 1 μ g, Lane 3: 2 μ g, Lane 4: 5 μ g Cyclophilin A.

Handling & Storage

Use/Stability	Caution: Loss of protein may occur at high dilutions due to adsorption to container surfaces. Freezing of the enzyme or storage of the enzyme in diluted form could lead to loss of activity. In order to best maintain the endotoxin specification and peptidyl prolyl isomerase activity, it is recommended that aliquots be withdrawn with sterile pipet tips, if possible under a sterile hood.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Activity	Catalyzes the cis-trans-isomeration of X-Pro-peptide bonds. The protein folding that is limited by the isomeration of X-Pro-peptide bonds can be accelerated by cyclophilin A. Also binds to cyclosporins and the cyclophilin A/cyclosporin A (Prod. No. BML-SE163).
Alternative Name	Peptidylproline cis-trans-isomerase
Application Notes	Catalyzes the <i>cis-trans</i> -isomerization of X-Pro-peptide bonds. Protein folding reactions that are limited by the isomerization of X-Pro-peptide bonds are accelerated by cyclophilin A. Cyclophilin A binds tightly to the immunosuppressive agent Cyclosporin A. This complex binds calcineurin (PP2B) and inhibits its phosphatase activity.
Endotoxin Content	Packaged under sterile conditions with a maximum endotoxin content of 0.1 EU/μg determined using the chromogenic Limulus amebocyte lysate (LAL) test.
Formulation	Liquid. In 20 mM TRIS-HCl, pH 7.8.
MW	18kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> .

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)