Cyclophilin A (human), (recombinant)

Ordering Information

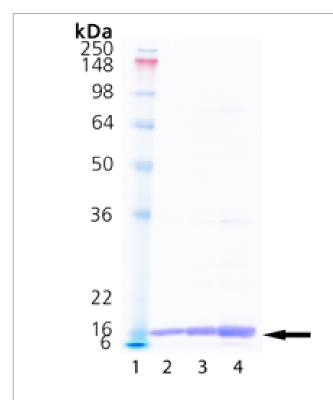
Order Online »

BML-SE105-0100

100µg

Manuals, SDS & CofA

View Online »



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

Handling & Storage

Use/StabilityCaution: 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to best maintain the endotoxin specification and peptidyl prolyl isomerase activity, it is recommended that aliquots be withdrawn with sterile pipet tips, if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 *cis-trans*-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 E. coli.

UniProt ID P62937

Last modified: May 29, 2024

