

# TCP-1 $\alpha$ monoclonal antibody (23c)

TCP-1 $\alpha$  is a 60 kDa subunit of a cytosolic hetero-oligomer chaperone that is known to be involved in the folding of Actin and Tubulin. This protein is a member of the Chaperonin family, which includes *E. coli* GroEL, the mitochondrial Heat-shock protein Hsp60, the plastid Rubisco-subunit-binding protein, and the archaeobacterial protein TF55. These Chaperonins assist the folding of proteins upon ATP hydrolysis.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 7

[View Online »](#)

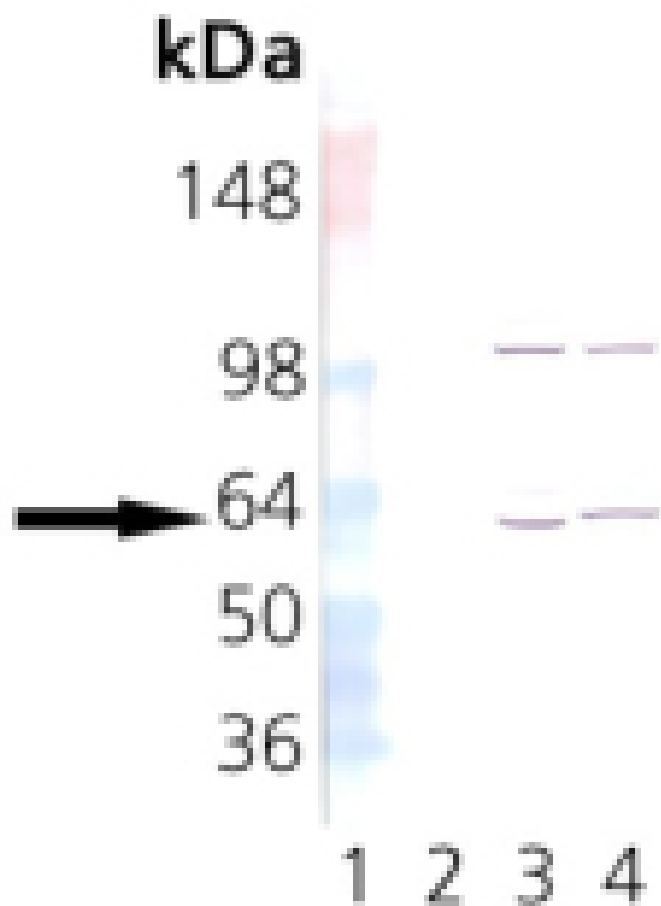
## Ordering Information

[Order Online »](#)

ADI-CTA-123-D	50 $\mu$ g
ADI-CTA-123-F	200 $\mu$ g

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of TCP-1a: Lane 1: MW marker, Lane 2: Hsp60 Recombinant Human Protein, Lane 3: 3T3 Heat Shocked, Lane 4: PC-12 Heat Shocked.

# Handling & Storage

Handling                      Avoid freeze/thaw cycles.

Long Term Storage           -20°C

Shipping                      Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name    CCT, p63, Tric, T-complex 1

Application    ICC, IHC, IP, WB

Application Notes    Detects a band of ~60kDa by Western blot.

Clone    23c

Formulation    Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

GenBank ID    D90344

Host    Rat

Immunogen    Recombinant mouse TCP-1α (C-terminal fragment).

Isotype    IgG2c

Purity Detail    Protein G affinity purified.

Recommendation Dilutions/Conditions    Western Blot (1:1,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source    Purified from mouse ascites.

Species Reactivity    Chicken, Dog, Hamster, Human, Mouse, Plant, Rabbit, Rat, Sheep

UniProt ID

P11983

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

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)