

TCP-1 α monoclonal antibody (23c)

TCP-1 α 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 »](#)

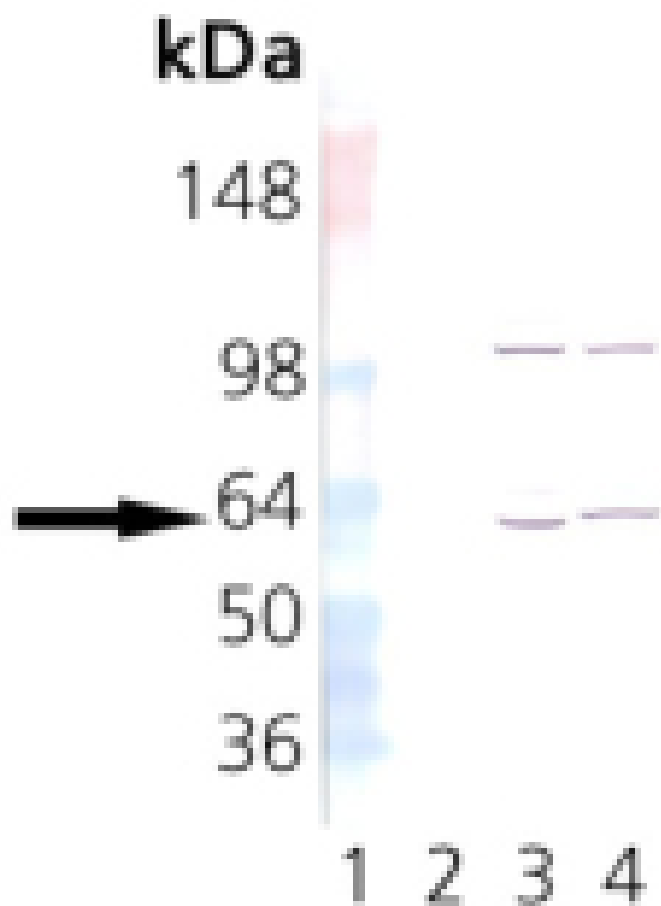
Ordering Information

[Order Online »](#)

ADI-CTA-123-D	50 μ g
ADI-CTA-123-F	200 μ 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 Western Blot (1:1,000, colorimetric) Suggested dilutions/conditions may not be available
Dilutions/Conditions 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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

P11986
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-
fr@enzolifesciences.com](mailto:info-fr@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)

Worry-free Guarantee

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

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.