

TCP-1 α monoclonal antibody (91a)

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: 3

[View Online »](#)

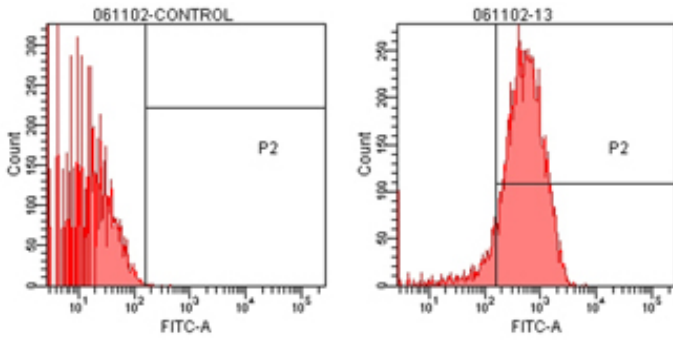
Ordering Information

[Order Online »](#)

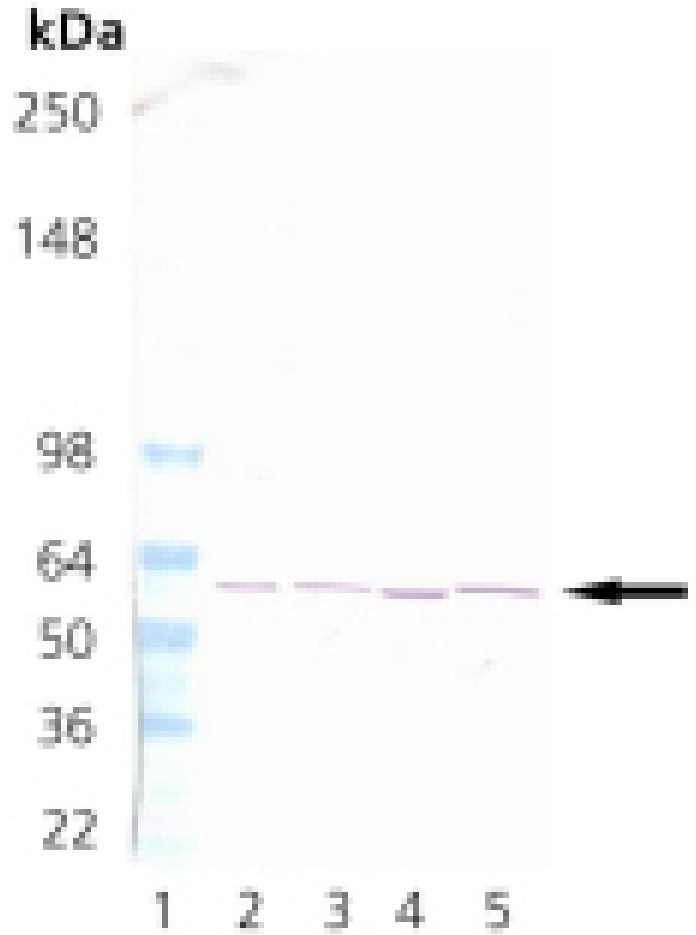
ADI-CTA-191-D	50 μ g
---------------	------------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis human colon cancer Coca-2 cells were analyzed by FLOW cytometry using isotype control antibody (left) or TCP-1alpha antibody, clone #91a (right).



Western blot analysis of CTA-191: Lane 1: MW marker, Lane 2: HeLa Cell Lysate, Heat shocked, Lane 3: PC-12 Cell Lysate Heat Shocked, Lane 4: 3T3 Cell Lysate, Heat Shocked, Lane 5, RK-13 Cell Lysate, 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 Flow Cytometry, ICC, IHC, IP, WB

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

Clone 91a

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

GenBank ID D90344

Host Rat

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

Isotype IgG2a

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:100)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 Bovine, C. elegans, Dog, Drosophila, Guinea pig, Hamster, Human, Monkey, Mouse, Plant, Porcine, Rabbit, Rat, Yeast

Technical Info / Product Notes

Recommended by the Human Protein Atlas Organization for IHC (Ensembl No. ENSG00000120438).

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)