

14-3-3η polyclonal antibody

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

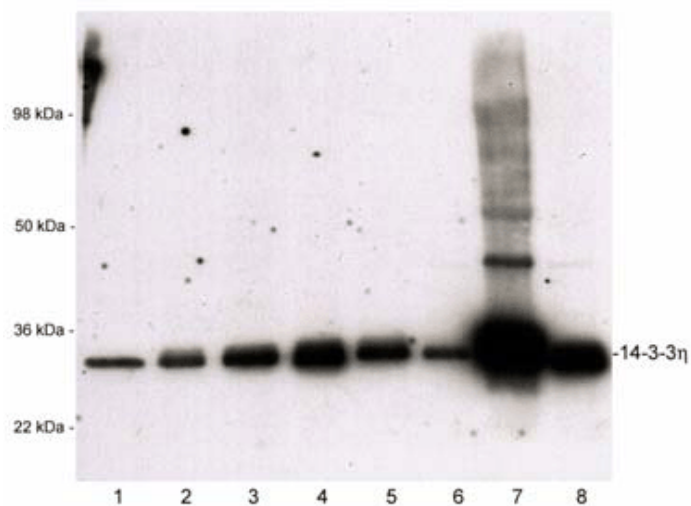
Ordering Information

[Order Online »](#)

BML-SA476-0100	100μl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot of 14-3-3 eta in 1) HeLa S100 (C), 2) HeLa whole cell, 3) Jurkat, 4) MDCK, 5) A431, 6) MDF-7, 7) rat brain, and 8) PC12 cell lysates. Developed with anti-14-3-3 eta antibody BML-SA476 (1:3'000) and HRP/ECL.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name KCIP-1, Protein kinase C inhibitor protein 1, Protein AS1

Application IHC, WB

Formulation Liquid. Neat serum containing sodium azide.

Host Rabbit

Immunogen Acetylated N-terminal domain of 14-3-3 eta.

Recommendation Dilutions/Conditions Immunohistochemistry (1:600)Western Blot (1:3,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Bovine, Chicken, Human, Mouse, Rat, Sheep

UniProt ID Q04917 (human)

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-be@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)