

14-3-3 ϵ (CT)

polyclonal antibody

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

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SA475-0100	100 μ l
----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

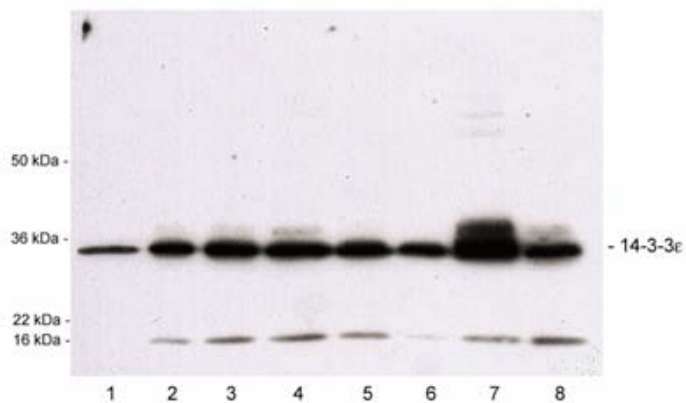


Figure: Western blot analysis of cell lysates using 14-3-3 ε, pAb (Prod. No. BML-SA475). Antibody dilution 1:3'000, blot was developed by HRP/ECL. 1. HeLa S100 (cytosol) 2. HeLa (whole cell) 3. Jurkat 4. MDCK 5. A4316. MDF-77. rat brain 8. PC12

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 KCIP-1, Protein kinase C inhibitor protein 1, 14-3-3E

Application ELISA, IHC, WB

Formulation Liquid. Neat serum. Contains no preservatives.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 239-255 (CG239EEQNKEALQDVEDENQ255) of 14-3-3ε.

Recommendation Western Blot (HRP/ECL detection, 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, Rabbit, Rat, Sheep

Specificity Specific for the ε isoform.

UniProt ID P62258 (human)

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



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)