

[pSer185]14-3-3 polyclonal antibody

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

Citations: 1

[View Online »](#)

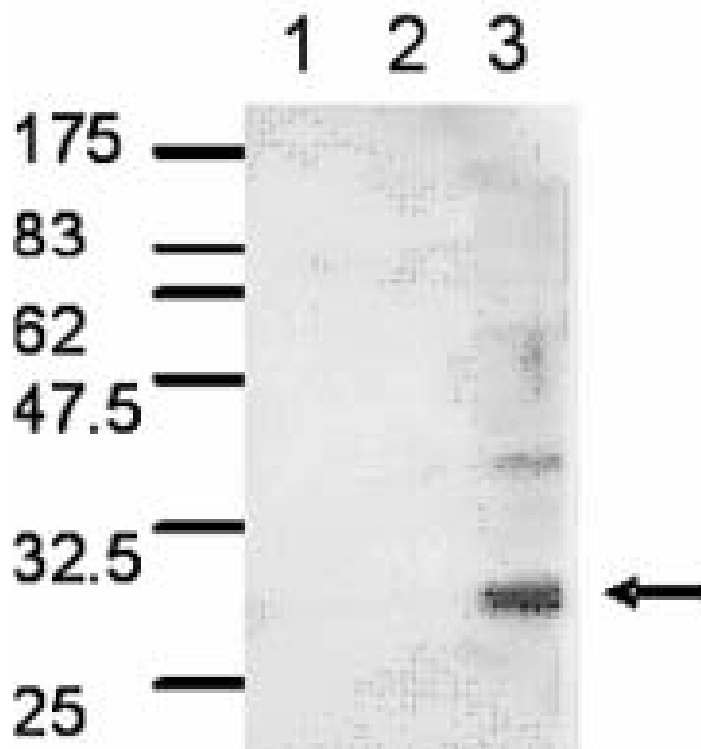
Ordering Information

[Order Online »](#)

BML-SA479-0025	25µl
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis performed with BML-SA479 antibody (1:500), secondary anti-sheep antibody (1:2'000), blotted against 2μg each of: 1. Recombinant 14-3-3β expressed in E.coli 2. Recombinant 14-3-3ζ expressed in E.coli 3. Sheep brain homogenate, which contains ~50% of β and ζ isoforms that are phosphorylated on Ser¹⁸⁵.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name KCIP-1, Protein kinase C inhibitor protein 1, 14-3-3 β / ζ , Protein 1054

Application WB

Host Sheep

Immunogen 14-3-3 phosphorylated at Ser185.

Purity Detail Affinity purified phospho-specific antibody.

Recommendation Western Blot (1:300-500) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Bovine, Human, Mouse, Rat, Sheep

Specificity Recognizes 14-3-3 β and 14-3-3 ζ phosphorylated at Ser¹⁸⁵.

UniProt ID P31946 (human 14-3-3 β), P63104 (human 14-3-3 ζ)

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)