

# [pSer185]14-3-3 polyclonal antibody

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

Citations: 1

[View Online »](#)

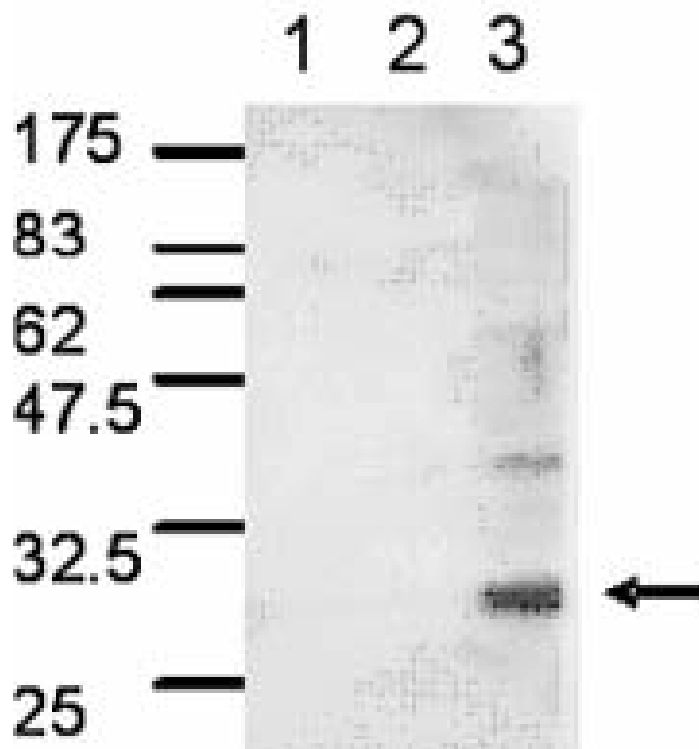
## Ordering Information

[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 Ser<sup>185</sup>.

## 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, 14-3-3β/ζ, Protein 1054
Application	WB
Host	Sheep
Immunogen	14-3-3 phosphorylated at Ser185.
Purity Detail	Affinity purified phospho-specific antibody.
Recommendation Dilutions/Conditions	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 14-3-3β and 14-3-3ζ phosphorylated at Ser <sup>185</sup> . .
UniProt ID	P31946 (human 14-3-3β), P63104 (human 14-3-3ζ)
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .

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-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)