

Dishevelled-2 polyclonal antibody

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

Citations: 1

[View Online »](#)

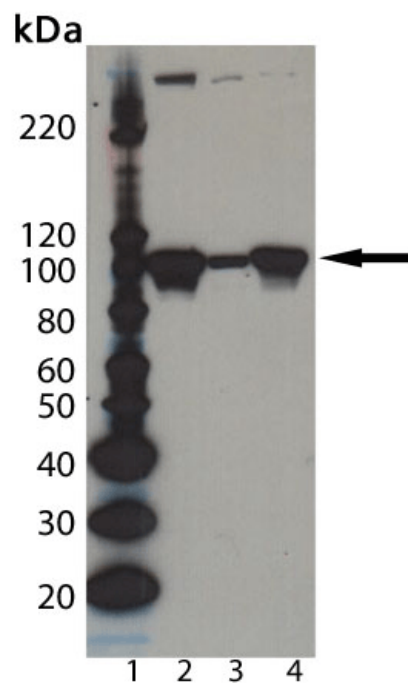
Ordering Information

[Order Online »](#)

BML-DA4270-0025	25µl
BML-DA4270-0100	100µl

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of Dishevelled-2 pAb (Prod. No. BML-DA4270): Lane 1: MW Marker, Lane 2: PC-3 cell lysate, Lane 3: Jurkat cell lysate, Lane 4: PC-12 cell lysate.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	DVL-2
Application	IHC, WB
Formulation	Liquid. In PBS, pH 7.4, containing 10mM sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 600-614 (CG600AGRTGRPEERAPES614) of human and mouse dishevelled 2 protein. The amino acid sequence does not exist in human or mouse dvl-1 or dvl-3.
Source	Purified from rabbit serum.
Species Reactivity	Human, Mouse
UniProt ID	O14641 (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-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)