

Blocking peptide for Dishevelled-2 pAb (Prod. No. BML- DA4270)

Sequence derived from human/mouse dishevelled-2 [Cys+600-614]. Used to generate the rabbit polyclonal antiserum to dishevelled-2 (BML-DA4270). The peptide is suitable for use in negative control (antibody blocking) experiments in Western blotting and immunohistochemistry.

Ordering Information

[Order Online »](#)

BML-DP9427-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance White powder.

MW 1671.8

Purity ≥86% (HPLC)

Sequence H-Cys-Gly-Ala-Gly-Arg-Thr-Gly-Arg-Pro-Glu-Glu-Arg-Ala-Pro-Glu-Ser-OH

Solubility Soluble in water and aqueous buffers.

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)