## Dishevelled-2 polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

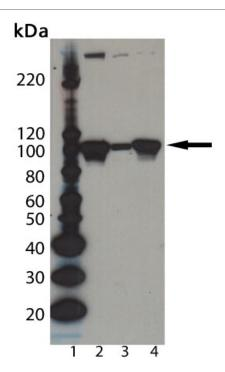
## **Ordering Information**

Order Online »

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

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

