

Biliverdin Reductase

polyclonal antibody

The phosphoprotein BVR (Biliverdin Reductase) is an oxireductase which catalyzes the reduction of the gamma-meso bridge of Biliverdin to Bilirubin. BVR is unique among all enzymes characterized to date because BVR has dual pH/dual cofactor requirements. The reductase is highly conserved in its primary structure and molecular properties and uses NADH in the acidic range (peak activity at pH~6.7), whereas NADPH is utilized in the basic range (peak activity at pH ~8.7).

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

Citations: 7

[View Online »](#)

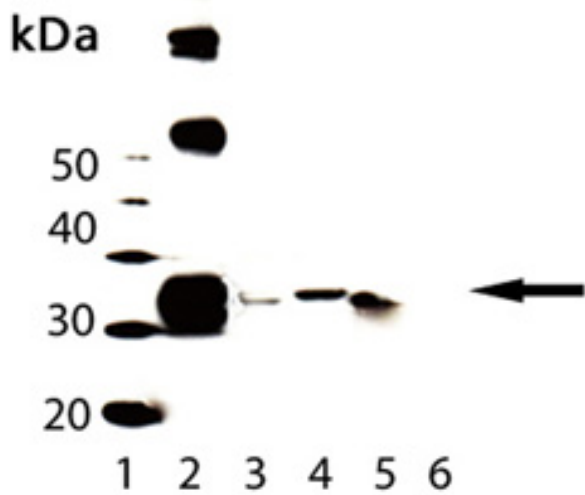
Ordering Information

[Order Online »](#)

ADI-OSA-450-E	100µl
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Biliverdin Reductase, pAb (Prod. No. ADI-OSA-450): Lane 1: MW Marker, Lane 2: Biliverdin Reductase Protein (rat) (recombinant) (Prod. No. ADI-OSP-450), Lane 3: Rat-2 Cell Lysate, Lane 4: 3T3 Cell Lysate (Prod. No. ADI-LYC-3T100), Lane 5: PC-12 Cell Lysate (Prod. No. ADI-LYC-PC100), Lane 6: HO-1 Protein (rat) (recombinant) (Prod. No. ADI-SPP-730)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name BLVR

Application IP, WB

Application Notes Detects bands of ~33kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

GenBank ID M81681

Host Rabbit

Immunogen Recombinant rat Biliverdin Reductase.

Purity Detail Ammonium sulfate precipitation purified.

Recommendation Dilutions/Conditions Immunoprecipitation (1:100)Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity E. coli, Mouse, Rat

UniProt ID P46844

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-
fr@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)

Last modified: May 29, 2024