

# Biliverdin Reductase (rat), (recombinant)

**Recombinant rat Biliverdin Reductase that catalyzes the reduction of biliverdin to bilirubin and modulates heme metabolism.**

Biliverdin reductase (BVR), Heme oxygenase (HO) and UDP-glucuronosyltransferases (UGTs) form the major erythrocytic heme catabolic pathway in humans and most mammalian species. BVR is an oxireductase which catalyzes the reduction of the gamma-meso bridge of biliverdin to bilirubin. Data shows BVR as a phosphoprotein, and establishes phosphorylation as essential to the activity of BVR to reduce biliverdin to bilirubin. These studies also found that the enzyme is autophosphorylated and that phosphorylation is reversible. In mammals, bilirubin acts as either a neurotoxin or an antioxidant depending on its ratio to protein and concentration, possesses immunomodulatory properties, induces Arylhydrocarbon receptor-mediated activation of cytochrome P450 1a1 expression, and inhibits protein phosphorylation. 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, using 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). A single copy gene encodes BVR, and expression of the ~1.6kb BVR message occurs abundantly in kidney, spleen, liver and brain, as well as at lower levels in the thymus and minimal levels in testis. The human BVR was found to resolve into four isoelectric zones and two molecular mass forms (40.7kDa and 39.6kDa) in the liver, whereas researchers detected variants in the kidney.

Citations: 1

[View Online »](#)

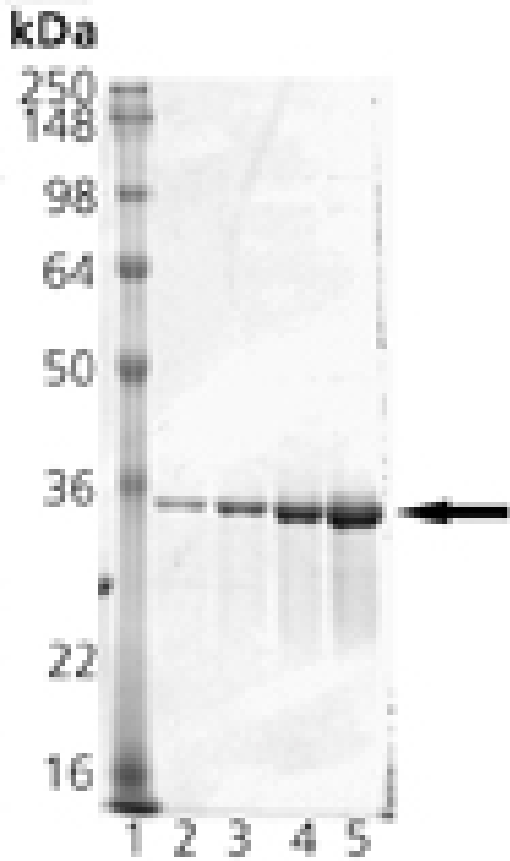
## Ordering Information

[Order Online »](#)

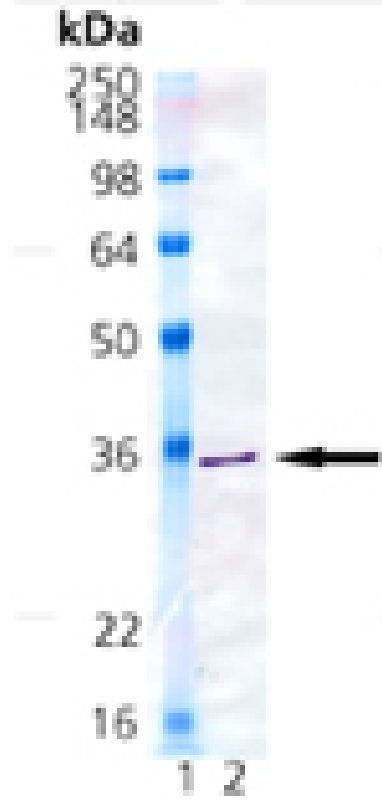
ADI-OSP-450-D	50µg
---------------	------

## Manuals, SDS & CofA

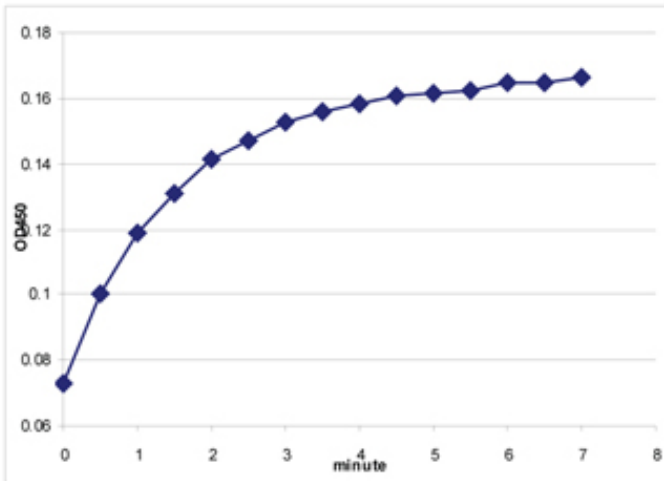
[View Online »](#)



SDS-PAGE analysis: Lane 1: MWM, Lane 2: 0.5 $\mu$ g, Lane 3: 1.0 $\mu$ g, Lane 4: 2.0 $\mu$ g, Lane 5: 5.0 $\mu$ g of Biliverdin Reductase Active Recombinant Protein.



Western Blot Analysis: Lane 1: MWM, Lane 2: Biliverdin Reductase Active Recombinant protein probed with Biliverdin Reductase pAb (Prod. No. ADI-OSA-400).



Biliverdin Reductase Activity: 1 $\mu$ g of Biliverdin Reductase Active Recombinant Protein was incubated with biliverdin in the presence of NADPH at pH 8.5, 37C. Product formation was monitored with a spectrometer at 450 nm.

## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	BLVR
Application Notes	Biliverdin Reductase Activity assay (positive). Western blot control.
Formulation	Liquid. In 50mM TRIS-HCl, pH 8.0, containing 1.0mM EDTA, and 5.0mM DTT.
GenBank ID	M81681
MW	~33kDa
Purity	≥95% (SDS-PAGE; Western blot)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> .
UniProt ID	P46844

Last modified: October 9, 2025



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)