

# LBP (mouse) monoclonal antibody (biG 33)

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

Citations: 4

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-804-502-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Long Term Storage            -20°C

Shipping                        Blue Ice

# Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Lipopolysaccharide binding protein
Application	ELISA, FUNC, WB
Application Notes	Functional Application: inhibits binding of LPS to membrane bound CD14.
Clone	biG 33
Crossreactivity	Cross-reacts weakly with rat LBP.
Formulation	Lyophilized. Contains PBS.
Host	Mouse
Immunogen	Recombinant mouse LBP (lipopolysaccharide binding protein) (in LBP knockout mouse).
Isotype	IgG1
Purity Detail	Protein-G affinity purified.
Recommendation Dilutions/Conditions	ELISA, Western Blot, Functional Application (Inhibits binding of LPS to membrane bound CD14)
Reconstitution	Reconstitute with 40µl distilled water.
Species Reactivity	Mouse
Technical Info / Product Notes	Binding titer at mouse LBP-ELISA: 1:50'000. Inhibits LPS-binding (0.5µg/ml) to CD14 <sup>+</sup> CHO cells (2×10 <sup>6</sup> /ml) up to 1:1'000, total inhibition is 1:5.

UniProt ID

Q61805

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)