

# LBP (mouse) monoclonal antibody (biG 33)

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

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-5020	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Lipopolysaccharide binding protein
<b>Application</b>	ELISA, FUNC, WB
<b>Application Notes</b>	Functional Application: inhibits binding of LPS to membrane bound CD14.
<b>Clone</b>	biG 33
<b>Crossreactivity</b>	Cross-reacts weakly with rat LBP.
<b>Formulation</b>	Lyophilized. Contains PBS.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant mouse LBP (lipopolysaccharide binding protein) (in LBP knockout mouse).
<b>Isotype</b>	IgG1
<b>Purity Detail</b>	Protein-G affinity purified.
<b>Recommendation Dilutions/Conditions</b>	ELISA, Western Blot, Functional Application (Inhibits binding of LPS to membrane bound CD14)
<b>Reconstitution</b>	Reconstitute with 40µl distilled water.
<b>Species Reactivity</b>	Mouse
<b>Technical Info / Product Notes</b>	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)