

CD14 (mouse) monoclonal antibody (biG 53)

The CD14 glycoprotein, gp 55, is present on most monocytic and macrophage like cell types: monocytes, macrophages, Kupffer cells, pleural phagocytic cells and dendritic reticular cells. CD14 is also observed on granulocytes and activated or transformed B-cells. Furthermore CD14 is present in a soluble form in mouse serum, urine and other body fluids. The CD14 molecule has been reported to be a receptor for endotoxin.

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

Ordering Information

[Order Online »](#)

ALX-804-499-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, Flow Cytometry, FUNC

Application Notes Functional Application: inhibits binding of LPS to CD14; up to 1µg/ml CD14 inhibits binding of 0.5µg/ml FITC-LPS to 6×10^5 CD14⁺ CHO transfectants (FACS).

Clone biG 53

Formulation Lyophilized. Contains no preservatives.

Host Mouse

Immunogen Recombinant mouse CD14.

Isotype IgG1

Purity Detail Protein G-affinity purified.

Reconstitution Reconstitute with 100µl distilled water.

Species Reactivity Mouse

Specificity Binds to LPS-activated mouse monocytes.

Technical Info / Product Notes Binding titer at mouse CD14 transfected CHO-cells is more than 20'000.

UniProt ID P10810

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)