

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](#)

.



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)