

CD14 monoclonal antibody (biG 10)

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 human 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](#).

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-496-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, WB

Application Notes Functional Application: Inhibits binding of LPS to CD14; total inhibition at 100µg/ml.

Clone biG 10

Formulation Lyophilized from PBS. Contains no preservatives.

Host Mouse

Immunogen Monocytes and immunoaffinity purified soluble human CD14.

Isotype IgG1

Purity Detail Protein G affinity-purified.

Reconstitution Reconstitute with 100µl distilled water.

Species Reactivity Bovine, Dog, Horse, Human, Porcine

Specificity Recognized epitopes in human: aa 9-13 and aa 39-44.

Technical Info / Product Notes Binding titer at human CD14 transfected CHO-cells is more than 10'000.

UniProt ID P08571

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)