

CD14, soluble (human) ELISA kit

The CD14, soluble (human) ELISA kit is a complete assay for the quantitative determination of soluble human CD14 concentration. This kit is to be used with serum, plasma and cell culture medium from human origin.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-302-KI01	96 wells
------------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	For long term storage keep vials 3 and 4 at -20°C. Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ELISA
Application Notes	For the quantitative determination of soluble human CD14 in serum, plasma and cell culture medium.
Assay Time	2 hours + 13+/- 1 minute
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	1 precoated ELISA module 1 vial detecting antibody (POD-labelled monoclonal antibody to human CD14) (ready-to-use) 1 vial human CD14 standard (recombinant human CD14, lyophilized) 1 vial reference serum (lyophilized) 2 tablets PBS 1 vial dilution buffer 1 vial Tween 20 1 vial stopping solution (ready-to-use) 1 vial substrate solution (ready-to-use)
Quantity	96 wells (~80 tests)
Sensitivity	Range: 5-50ng/ml
Species Reactivity	Human
UniProt ID	P08571
Wavelength	450 nm



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)