

SDMA ELISA kit

The SDMA ELISA kit is a complete assay for the quantitative determination of symmetric dimethylarginine (SDMA) levels. This kit is to be used with EDTA-plasma and serum from human origin. The SDMA ELISA has no cross-reactivity with ADMA, NMMA or L-arginine.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-331-KI01	96 wells
------------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Store standard and controls at -20°C. Store rest of the kit at +4°C.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NG, N'-G-Dimethyl-L-arginine ELISA kit

Application ELISA

Application Notes For the quantitative determination of symmetric dimethylarginine (SDMA) in human EDTA-plasma and serum. Does not cross-react with ADMA, NMMA or L-arginine.

Assay Time 15 to 20 hours + 1 hour

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Quantity 96 wells (~80 tests).

Sensitivity 0.05µmol/l (range 0.1 to 2µmol/l, expected value: 0.47µmol/l ± 0.14µmol/l)

Species Reactivity Human

Wavelength 450 nm

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-
fr@enzolifesciences.com](mailto:info-fr@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)