

ADMA direct (mouse/rat) ELISA kit

The ADMA direct (mouse/rat) ELISA kit is a complete assay for the quantitative determination of ADMA concentration. This kit is to be used with EDTA-plasma, serum, and cell culture supernatants from mouse or rat origins. The ADMA direct (mouse/rat) ELISA has no cross-reactivity with SDMA, NMMA or L-arginine.

Asymmetric dimethylarginine (ADMA) is an endogenous inhibitor of nitric oxide (NO) synthesis. NO is one of the major endothelium-derived vasoactive mediators and is involved in the modulation of blood flow and blood pressure. Even small changes of ADMA concentration alter vascular NO production, vascular tone, and systemic vascular resistance. Elevated ADMA concentrations are found in patients with diabetes mellitus, hypercholesterolemia, hypertension, pulmonary hypertension, peripheral arterial disease, chronic renal failure, preeclampsia or erectile dysfunction. ADMA is therefore a cardiovascular risk factor.

Citations: 17

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-327-KI01	96 wells
------------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA

Application Notes For the quantitative detection of asymmetrical dimethylarginine (ADMA) in mouse or rat EDTA-plasma, serum, and cell culture supernatants. Does not cross-react with SDMA, 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

Species Reactivity Mouse, Rat

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)