

# ADMA (human) ELISA kit

The ADMA (human) ELISA kit is a complete assay for the quantitative determination of ADMA concentration. This kit is to be used with serum, citrate- and EDTA-plasma from human origin. The ADMA (human) 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: 11

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-850-323-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 determination of asymmetrical dimethylarginine (ADMA) in human serum, citrate- and EDTA-plasma. 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 (expected value: 0.45µmol/l ± 0.19µmol/l)

**Species Reactivity** Human

**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)