

# 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

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-  
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)