

# 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 »](#)

