

Endothelin-1 ELISA kit

Sensitive, rapid Endothelin-1 ELISA kit for cardiovascular and stress-related research.

The Endothelin-1 EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 2 hours. This ultra-sensitive kit lets you quantify as little as 0.41 pg/mL ET-1 in a variety of biological matrices with the option of overnight or a 1 hour primary incubation.

Endothelins (ET) are isopeptides produced by vascular endothelium that have potent vasoconstrictive activity. The peptides are encoded by three separate genes and processed to yield 39-residue "Big ET" molecules which are further processed to the 21 amino acid sequences designated ET-1, ET-2 and ET-3. ET action is mediated via binding to two receptors, ET_A and ET_B. Activation of ET_A increases vasoconstriction and sodium retention, thereby increasing blood pressure. Activation of ET_B receptors triggers the release of vasorelaxive factors such as nitric oxide (NO) and prostanooids from endothelial cells and increases urine production and elimination, thereby lowering blood pressure.

Citations: 60

[View Online »](#)

Ordering Information

[Order Online »](#)

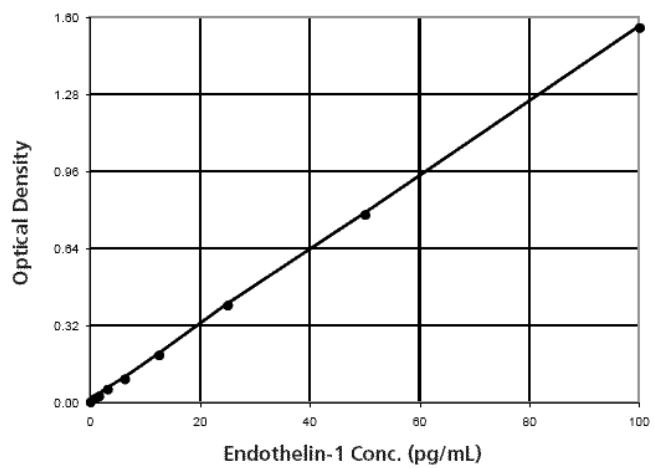
ADI-900-020A

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive measurement, quantify as little as 0.41 pg/ml ET-1, allowing you read both normal and pathological samples
- Higher throughput format with results in just 2 hours for up to 38 samples in duplicate
- Flexible option of overnight or 1 hour primary incubation
- Suitable for a variety of sample matrices



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of ET-1 in culture supernatants, plasma, and serum of human, mouse, rat, cow, dog, pig, and rabbit origin. Cited sample types include cell lysates and tissue homogenates.

Assay Time 2 hours

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

Contents Microtiter plate, Antibody diluent, Antibody concentrate, Assay buffer 17, Wash buffer concentrate, Standard, TMB Substrate, Stop solution 2

Sensitivity 0.41 pg/ml (range 0.78 – 100 pg/ml)

Species Reactivity Bovine, Dog, Human, Mouse, Porcine, Rabbit, Rat

UniProt ID P05305

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-
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)