

# Dopamine ELISA Kit

**Sensitive ELISA kit for the quantification of Dopamine in any species**

The Dopamine ELISA kit is a colorimetric immunoassay kit with results in < 2 hours. Absorbance is read at 450 nm. Measure up to 40 samples in duplicate per 96 well kit.

**Citations:** 9

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-KIT188-0001	96 wells
-----------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- High-sensitivity ELISA enables quantification of low amounts of Dopamine
- Rapid quantitative results in <2 hours
- High-throughput immunoassay measures up to 40 samples in duplicate per 96 well kit
- Pre-coated plates reduce errors and save time

## 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 dopamine in serum, plasma, tissue homogenates, and other biological fluids.
Assay Time	< 2 hours
Compatibility	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
Contents	Microtiter plate, Lyophilized Standard, Sample/standard dilution buffer, Concentrated biotin detector antibody, Antibody Dilution Buffer, HRP-Streptavidin Conjugate, Conjugate Dilution Buffer, TMB Substrate, Stop Solution, Wash Buffer (25X), Plate Sealer
Crossreactivity	Negligible crossreactivity to similar biological molecules was observed.
Sensitivity	0.938 ng/mL (1.56 – 100 ng/mL)
Species Reactivity	Species independent
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-  
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)