

NGAL (dog) ELISA kit

A colorimetric immunometric ELISA kit for the rapid measurement of NGAL in dog samples. Absorbance is read at 450nm.

NGAL (also known as Lipocalin-2 or Neutrophil Gelatinase-associated Lipocalin) is a small, ~25kDa secreted protein expressed in epithelial tissues throughout the body.

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

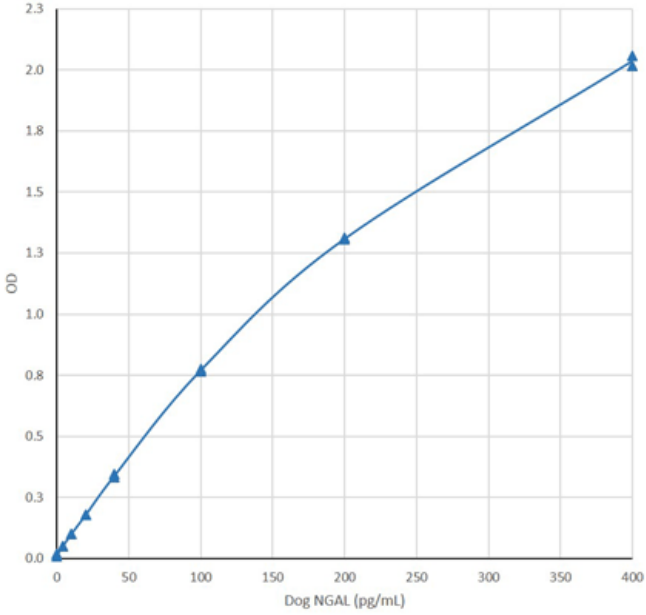
BPD-KIT-043	96 wells
-------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Ready-to-use - pre-coated plate and individual calibrators save time and reduce errors
- Fast - results from up to 40 samples in duplicate in just 4 hours per 96 well plate

Dog NGAL ELISA Kit (KIT 043) - Calibration curve



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Neutrophil gelatinase-associated lipocalin, Lipocalin 2, Siderocalin, Oncogene 24p3
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of NGAL in dog urine, plasma, serum, tissue extracts, or culture supernatants.
Assay Time	4 hours
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	1: Microwell plate coated with anti-NGAL (dog) antibody (96 wells) 2: 5X Sample Diluent 3: NGAL Dog Calibrators 1-8; 0, 4, 10, 20, 40, 100, 200, 400 pg/mL. 4: 25X Wash Solution Concentrate 5: Biotinylated Dog NGAL Antibody 6: HRP-Streptavidin 7: TMB Substrate 8: Stop Solution
Sensitivity	0.56pg/ml (range 4-400pg/ml)
Species Reactivity	Dog
Technical Info / Product Notes	For the manual please click here .
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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)