

# 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 »](#)

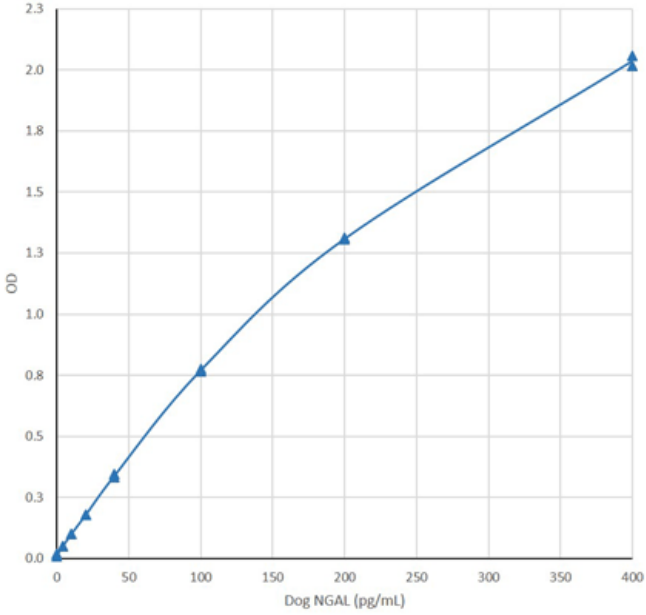
BPD-KIT-043	96 wells
-------------	----------

## 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



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)