

# NGAL (pig) ELISA kit

A colorimetric immunometric ELISA kit for the rapid measurement of NGAL in pig 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: 11

[View Online »](#)

## Ordering Information

[Order Online »](#)

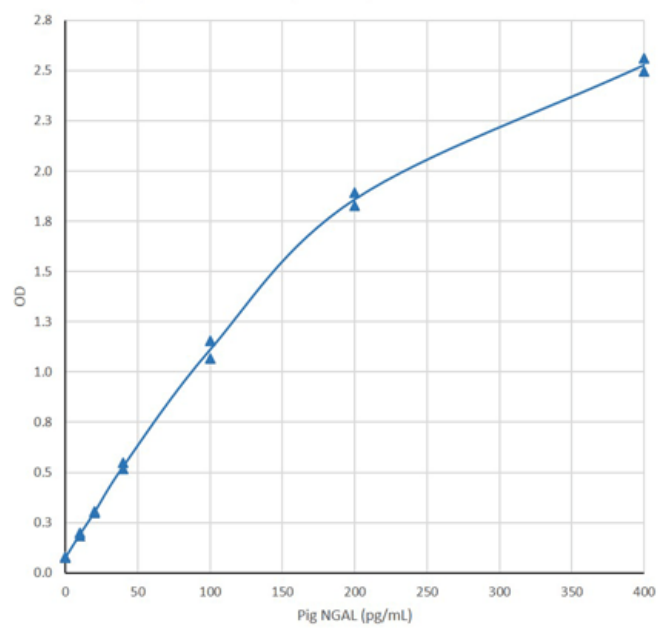
BPD-KIT-044	96 wells
-------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Ready-to-use - pre-coated plate and individual calibrators save time and reduce errors
- Easy-to-use - convenient storage at 4°C with room temperature incubations

Pig NGAL ELISA Kit (KIT 044) - 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 pig 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 (96 precoated wells); 2: 50ml 5x Sample Diluent Conc.; 3: 7 x 1ml Pig NGAL Calibrator (1-7. 0, 10, 20, 40, 100, 200, 400 pg/mL); 4: 40ml 25x Wash Solution Conc.; 5: 12ml Biotinylated Pig-NGAL Antibody; 6: 12ml HRP-Streptavidin; 7: 12ml TMB Substrate; and 8: 12ml Stop Solution.

**Sensitivity** 1pg/ml (range 10-400pg/ml)

**Species Reactivity** Porcine

**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)