

# NGAL (rat) ELISA kit

A colorimetric immunometric ELISA kit for the rapid measurement of NGAL in rat 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: 15

[View Online »](#)

## Ordering Information

[Order Online »](#)

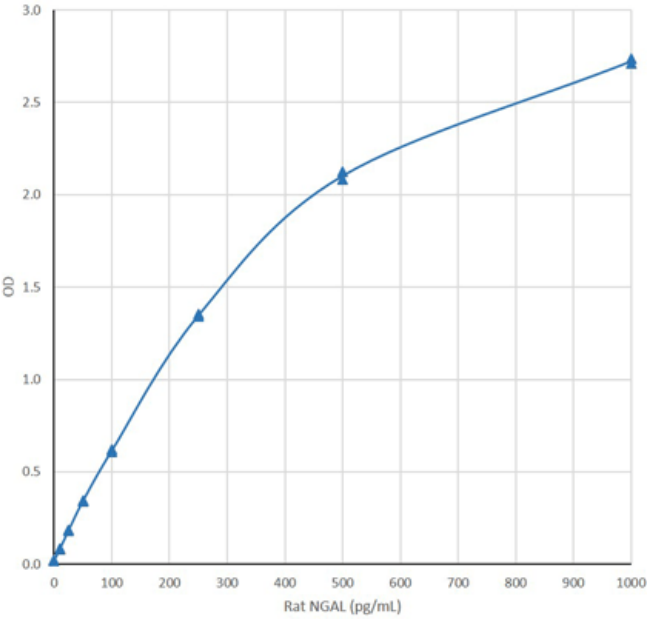
BPD-KIT-046	96 wells
-------------	----------

## 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

Rat NGAL ELISA Kit (KIT 046) - 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 rat urine, plasma, serum, tissue extracts, or culture supernatants.
Assay Time	4 hours
Compatibility	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
Contents	1: Microwell plate coated with anti-NGAL (rat) antibody (96 wells) 2: 5X Sample Diluent 3: NGAL Rat Calibrators 1-8. 4: 25X Wash Solution Concentrate 5: Biotinylated Rat NGAL Antibody 6: HRP-Streptavidin 7: TMB Substrate 8: Stop Solution
Sensitivity	0.5pg/ml (range 4-400pg/ml)
Species Reactivity	Rat
Technical Info / Product Notes	For the manual please <a href="#">click here</a> .
UniProt ID	P30152
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)