

# NGAL (mouse) ELISA kit

A colorimetric immunometric ELISA kit for the rapid measurement of NGAL in mouse 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: 8

[View Online »](#)

## Ordering Information

[Order Online »](#)

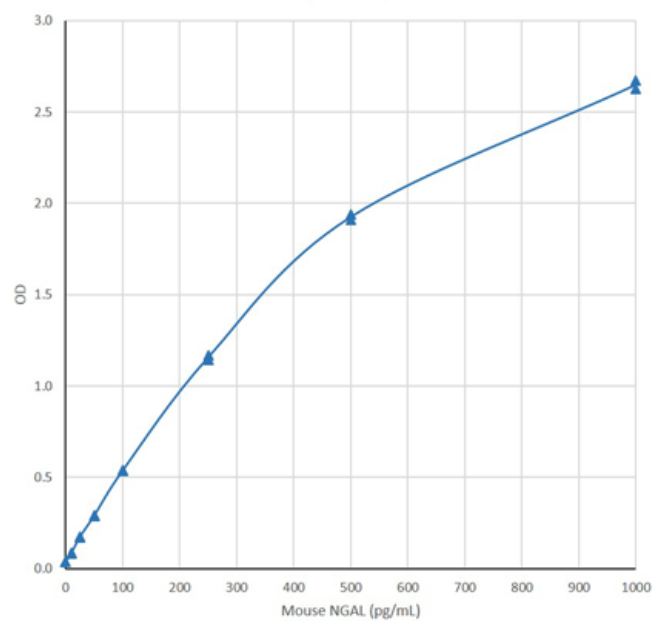
BPD-KIT-042	96 wells
-------------	----------

## 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

Mouse NGAL ELISA Kit (KIT 042) - 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 mouse 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 (mouse) antibody (96 wells) 2: 5X Sample Diluent 3: 2 containers of lyophilized Mouse NGAL calibrators. 4: 25X Wash Solution Concentrate 5: Biotinylated Mouse NGAL Antibody 6: HRP-Streptavidin 7: TMB Substrate 8: Stop Solution

### Sensitivity

0.75pg/ml (range 10-1000pg/ml)

### Species Reactivity

Mouse

### Technical Info / Product Notes

For the manual please [click here](#).

### UniProt ID

P11672

### Wavelength

450 nm



ENZO LIFE SCIENCES,  
INC.  
May 29, 2024  
info-  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
info-  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
info-  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
info-  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
info-  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
info-  
[uk@enzolifesciences.com](mailto:uk@enzolifesciences.com)