

# NGAL (human) ELISA kit

Highly sensitive ELISA, measuring as little as 4 pg/ml of NGAL

A colorimetric immunometric ELISA kit for the measurement of NGAL in human 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: 86

[View Online »](#)

## Ordering Information

[Order Online »](#)

BPD-KIT-036

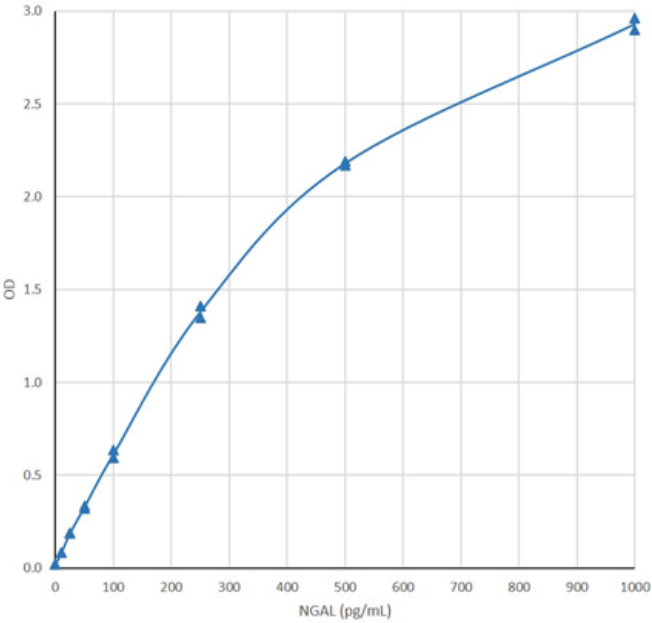
96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Fast results from up to 40 samples in duplicate in just 4 hours per 96 well plate
- Highly sensitive measurement, detecting as low as 4 pg/ml of NGAL
- Ready-to-use pre-coated plate and individual calibrators save time and reduce errors
- Fully quantitative results that surpass semi-quantitative Western blot analysis

Human NGAL ELISA Kit (KIT 036) - 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 human urine, plasma, serum, tissue extracts, and culture supernatants.

**Assay Time** <4 hours

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**Contents** 1: 12 strips of 8 Microwells each coated with anti-NGAL (96 wells) + frame 2: Sample Diluent 3: NGAL Calibrators 1-8; 0, 10, 25, 50, 100, 250, 500, 1000 pg/mL 4: 25X Wash Solution Concentrate 5: Biotinylated NGAL Antibody 6: HRP-Streptavidin Conjugate 7: TMB Substrate 8: Stop Solution

**Sensitivity** 4pg/ml (range 10-1000pg/ml)

**Species Reactivity** Human

**Technical Info / Product Notes** For the manual please [click here](#).

**UniProt ID** P80188

**Wavelength** 450 nm



Last modified: May 29, 2024

ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430

[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989

[info-be@enzolifesciences.com](mailto:info-be@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)