

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 »](#)

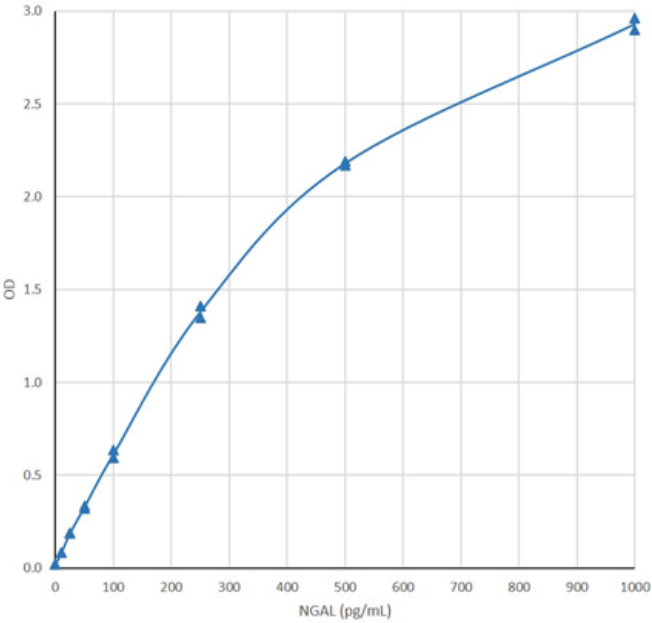
BPD-KIT-036	96 wells
-------------	----------

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



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)