

# NGAL (pig), (recombinant)

Pig neutrophil gelatinase-associated lipocalin (NGAL; also called lipocalin 2 or siderocalin) may, by analogy with its homologues in other mammalian species, be released from activated neutrophils in inflammation or infection, from certain epithelial cancers, and more dramatically, from renal tubular cells in response to ischemic or nephrotoxic injury.

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

Short Term Storage +4°C

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

**Concentration** 400pg/ml

**Formulation** Liquid. In 0.01 M phosphate buffer, pH 7.4, containing 0.14 M NaCl, 0.5% BSA, 0.5% Tween 20, 0.0016% Phenol red, and 0.1% Proclin 950.

**Source** Produced in *E. coli*.

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-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)