

NGAL (pig) monoclonal antibody (28)

Pig neutrophil gelatinase-associated lipocalin (NGAL; also 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.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

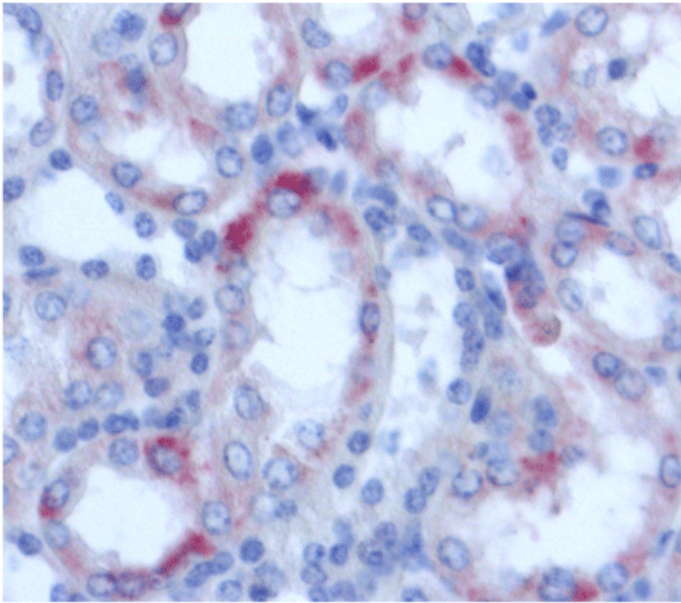
Ordering Information

[Order Online »](#)

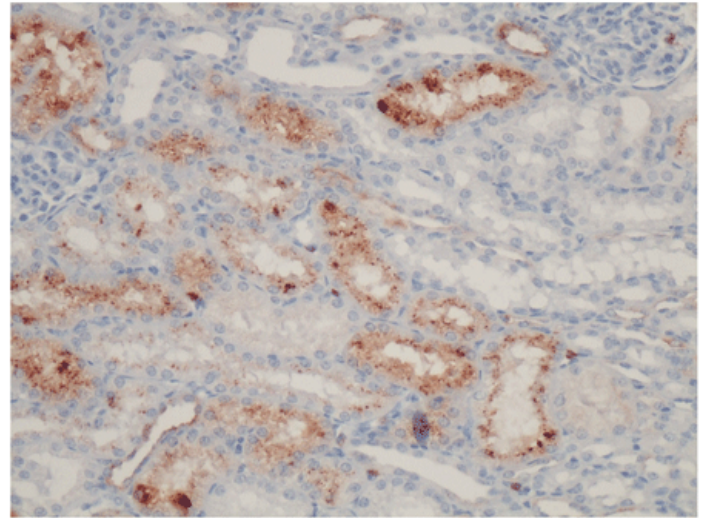
BPD-ABS-048-28-1	1mg
BPD-ABS-048-28-04	400µg

Manuals, SDS & CofA

[View Online »](#)



Positive staining of renal tubular cell cytoplasm, using mouse monoclonal antibody against pig NGAL.



NGAL expression in the ischemic injured pig kidney; detected with anti-pig NGAL antibody.

Handling & Storage

Short 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	ELISA, IHC (PS), WB
Application Notes	Detects a band of ~25kDa by Western blot.
Clone	28
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
Host	Mouse
Immunogen	Recombinant pig NGAL.
Isotype	IgG1κ
Purity Detail	Protein A/G-affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections 1:25)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Porcine
UniProt ID	F1RRX1
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)