

NGAL (dog) monoclonal antibody (17)

Dog 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](#).

Ordering Information

[Order Online »](#)

BPD-ABS-047-17-1	1mg
BPD-ABS-047-17-04	400µg

Manuals, SDS & CofA

[View Online »](#)

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, WB

Application Notes Detects a band of ~25kDa by Western blot.

Clone 17

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Recombinant dog NGAL.

Isotype IgG2ak

Purity Detail Protein A-affinity purified.

Species Reactivity Dog

UniProt ID E2RSM4

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)