

# NGAL (dog) monoclonal antibody (14)

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](#).

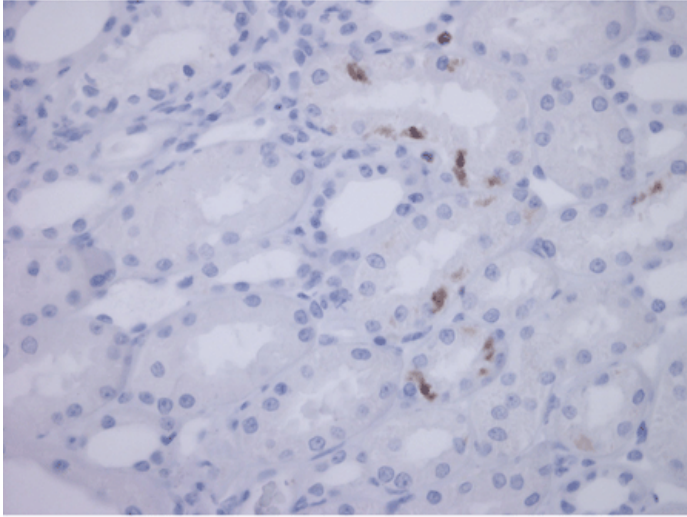
## Ordering Information

[Order Online »](#)

BPD-ABS-047-14-1	1mg
BPD-ABS-047-14-04	400µg

## Manuals, SDS & CofA

[View Online »](#)



IHC staining in dog kidney tissue. Scattered tubules cells were stained strongly from kidney with acute nephritis. Dilution: 1:50.

## 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, WB

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

**Clone** 14

**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** IgG2bk

**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)