

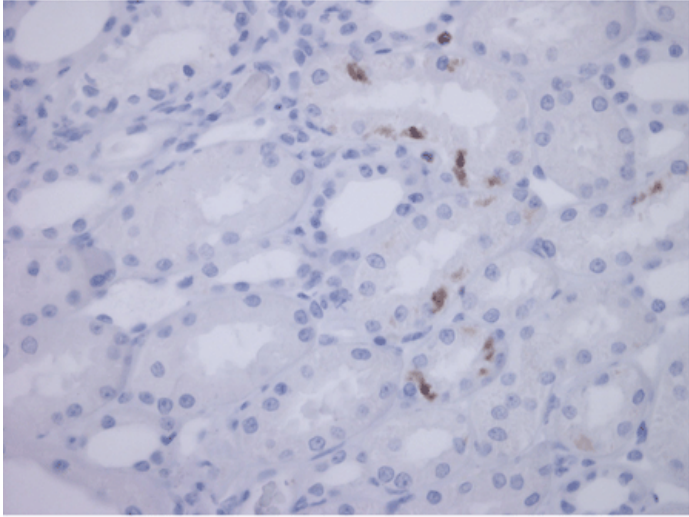
NGAL (dog) monoclonal antibody (14)

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](#).

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com