

# NGAL (mouse) monoclonal antibody (10)

Mouse neutrophil gelatinase-associated lipocalin (NGAL; also lipocalin 2, 24p3, SIP24, uterocalin and siderocalin) belongs to the lipocalin family of proteins which bind and transport small lipophilic molecules. Following research showing 24p3 dramatically upregulated in kidney cell cultures infected with SV40, it was subsequently identified as a lipocalin. Unlike human NGAL, mouse NGAL is an acute phase protein upregulated in the liver, but like its mammalian homologues, it is upregulated and secreted by the kidney in response to ischemic or nephrotoxic injury. It is also upregulated in the involuting uterus and mammary gland.

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

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	10
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
Host	Rat
Immunogen	Recombinant mouse NGAL.
Isotype	IgG1κ
Purity Detail	Protein A/G-affinity purified.
Species Reactivity	Mouse
UniProt ID	P11672
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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-fr@enzolifesciences.com](mailto:info-fr@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)