

NGAL (mouse) monoclonal antibody (29) (biotin conjugate)

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
Application Notes	Detects a band of ~25kDa by Western blot.
Clone	29
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.14mM sodium chloride and 15mM sodium azide.
Host	Rat
Immunogen	Recombinant mouse NGAL.
Isotype	IgG1k
Purity Detail	Protein A/G-affinity purified.
Species Reactivity	Mouse
UniProt ID	P11672
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