

NGAL (rat) monoclonal antibody (32) (biotin conjugate)

Rat lipocalin-2 is also called rat NGAL, as it is the ortholog of human neutrophil gelatinase-associated lipocalin (NGAL). Rat NGAL is a 25-kDa alpha 2-microglobulin-related protein or neu-related lipocalin. It is expressed by neutrophils and various epithelial cells, including cells of the renal proximal tubules. It is involved in physiological processes such as the binding of siderophore iron and tissue growth and regeneration. It is upregulated in a variety of pathological situations including infection, inflammation, ischemia-reperfusion and in certain adenocarcinomas.

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Neutrophil gelatinase-associated lipocalin, Lipocalin 2, Siderocalin, Neutrophil lipocalin

Application ELISA

Clone 32

Crossreactivity Does not cross-react with human, pig, dog or mouse.

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.14M sodium chloride and 15mM sodium azide.

GenBank ID 170496

Host Mouse

Immunogen Recombinant rat NGAL precursor.

Isotype IgG1κ

Purity Detail Protein A/G-purified.

Species Reactivity Rat

UniProt ID P30152

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



ENZO LIFE SCIENCES,
Last modified **May 29, 2024**
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