

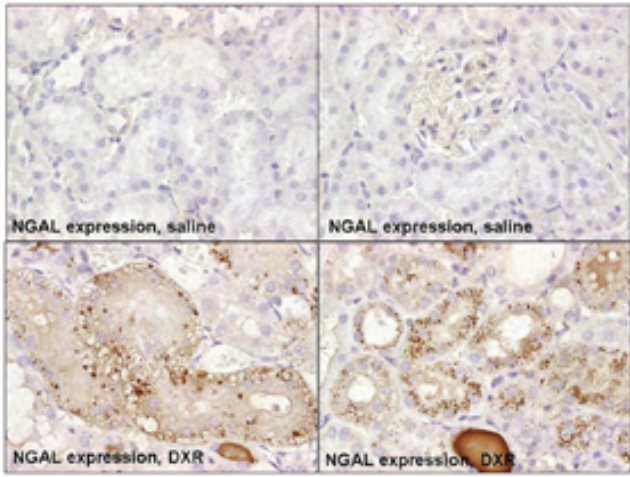
NGAL (rat) monoclonal antibody (8)

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 »](#)



Immunohistochemical analysis of kidney, formalin-fixed, paraffin-embedded tissue sections of cortical epithelial cells in the S1/S2 segments (proximal convoluted tubules) following DXR-induced acute kidney injury stained with NGAL (rat), mAb. Top panel, saline treatment; bottom panel, DXR treatment; X630.

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
Application Notes	Detects a band of ~25kDa by Western blot.
Clone	8
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
Host	Mouse
Immunogen	Recombinant rat NGAL precursor.
Isotype	IgG1κ
Purity Detail	Protein A/G-affinity purified.
Species Reactivity	Rat
UniProt ID	P30152
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-fr@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