

# 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](#).

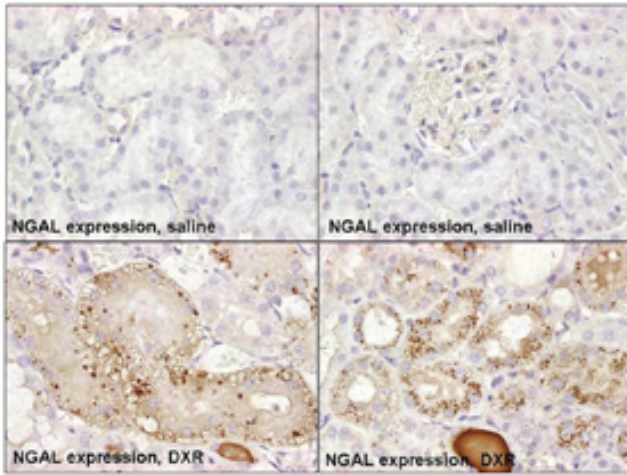
## Ordering Information

[Order Online »](#)

BPD-ABS-039-08-1	1mg
BPD-ABS-039-08-04	400µg

## 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](#).



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