

# NGAL (human) monoclonal antibody (14)

Neutrophil gelatinase-associated lipocalin (NGAL; also called lipocalin 2, siderocalin and neutrophil lipocalin) is a member of the lipocalin family of proteins which bind and transport small lipophilic molecules. NGAL is released by activated neutrophils and occurs as 25-kDa glycosylated single protein chain monomers, which form dimers and small amounts of higher oligomers, as well as complexes with matrix metalloproteinase 9 (MMP-9; gelatinase B). Low level expression of NGAL in a variety of epithelia may be increased in inflammation or cancers. Expression of NGAL in the kidney is dramatically increased by acute kidney injury.

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

## Ordering Information

[Order Online »](#)

BPD-ABS-062-14-1	1mg
BPD-ABS-062-14-04	400µg

## 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	14
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
GenBank ID	3934
Host	Mouse
Immunogen	Human recombinant NGAL coupled to carrier
Isotype	IgG2ak
Purity Detail	Protein A/G-purified.
Source	From culture supernatant.
Species Reactivity	Human
UniProt ID	P80188
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>