## NGAL (human) monoclonal antibody (15) (biotin conjugate)

Neutrophil gelatinase-associated lipocalin (NGAL; also called lipocalin 2, siderocalin and neutrophil lipocalin) belongs to 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-038-15B-015

150µg

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, WB

Clone 15

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

**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 Recombinant human NGAL.

**Isotype** IgG1κ

Purity Detail Protein A/G-purified.

Species Reactivity Human

UniProt ID P80188

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

