NGAL (human) monoclonal antibody (15)

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-038-15-1	1mg
BPD-ABS-038-15-04	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, 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 IgG1k

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 Worry-Free Guarantee



ENZO LIFE SCIENCES, Phone: 800.942.0430 info-

ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com

European Sales Office

eu@enzolifesciences.com

& Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Belgium, The Netherlands

France Phone: +33 472 440 655 fr@enzolifesciences.com

Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 uk@enzolifesciences.com