# NGAL (rat), (recombinant)

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.

### **Ordering Information**

Order Online »

ENZ-PRT115-1000

1ml

Manuals, SDS & CofA

View Online »

#### **Handling & Storage**

Short Term Storage +4°C

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, Oncogene 24p3

Concentration 400pg/ml

**Formulation** Liquid. In 0.01 M phosphate buffer, pH 7.4, containing 0.14

M NaCl, 0.5% BSA, 0.5% Tween 20, 0.0016% Phenol red,

and 0.1% Proclin 950.

**Source** Produced in *E. coli*.

UniProt ID P30152

Last modified: May 29, 2024



eu@enzolifesciences.com