

PRODUCT INFORMATION



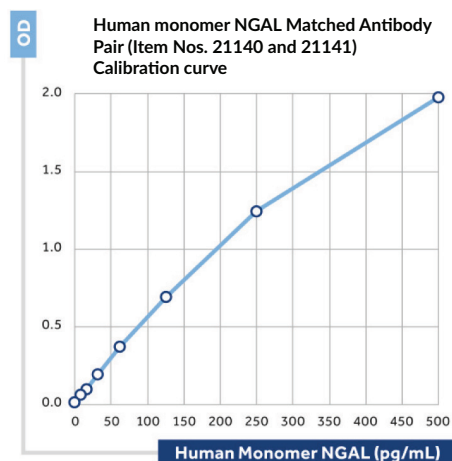
NGAL (human) Monoclonal Antibody (Clone 33)

Item No. 21141

Overview and Properties

Contents:	This vial contains 200 μ l or 1 ml of a 1 mg/ml (\pm 15%) solution of protein A-purified IgG.
Synonyms:	Lipocalin-2, Neutrophil gelatinase-associated lipocalin, Oncogene 24p3
Immunogen:	Human recombinant NGAL coupled to carrier and adsorbed to aluminum hydroxide gel
Species Reactivity:	(+) Human NGAL
Uniprot No.:	P80188
Form:	Liquid
Storage:	4°C (as supplied); store in the dark
Stability:	As supplied, 6 months from the QC date provided on the Certificate of Analysis, when stored properly
Storage Buffer:	0.01 M phosphate buffer, pH 7.4, containing 0.5 M sodium chloride and 15 mM sodium azide
Clone:	33
Host:	Mouse
Isotype:	IgG1/k
Application:	ELISA; NGAL (human) Monoclonal Antibody (Clone 33) binds human NGAL coated onto a solid phase. When it is used as a capture antibody with NGAL (human) Monoclonal Antibody - Biotinylated (Clone 14; Item No. 21140), the assay is specific for NGAL monomer, cross reacting <1% with NGAL homodimer.

Image



The figure illustrates the calibration curve of a sandwich assay with Item No. 21141 as the capture antibody and Item No. 21140 as the biotinylated detection antibody.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/22/2016

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Neutrophil gelatinase-associated lipocalin (NGAL) 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).¹ Low level expression of NGAL in a variety of epithelia may be increased in inflammation or cancers.² Expression of NGAL in the kidney is dramatically increased by acute kidney injury.

References

1. Kjeldsen, L., Johnsen, A.H., Sengeløv, H., *et al.* Isolation and primary structure of NGAL, a novel protein associated with human neutrophil gelatinase. *J. Biol. Chem.* **268**(14), 10425-10432 (1993).
2. Nielsen, B.S., Borregaard, N., Bundgaard, J.R., *et al.* Induction of NGAL synthesis in epithelial cells of human colorectal neoplasia and inflammatory bowel diseases. *Gut* **38**(3), 414-420 (1996).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM