

NKp46/NCR1

monoclonal antibody

(N1D9)

A natural cytotoxicity receptor (NCR), Nkp46, is a glycoprotein that has two extracellular Ig-like domains followed by a ~40 residue stalk region, a type I transmembrane domain, and a short cytoplasmic tail. Nkp46 has been shown to represent a novel NK cell-specific molecule involved in human NK cell activation. The natural cytotoxicity receptors (NCRs) are a recently characterized family of Ig-like activation receptors that appear to be major triggering receptors in tumor cell recognition. Nkp46 has been implicated in NK cell-mediated lysis of several autologous tumor cells and pathogen-infected cell lines.

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

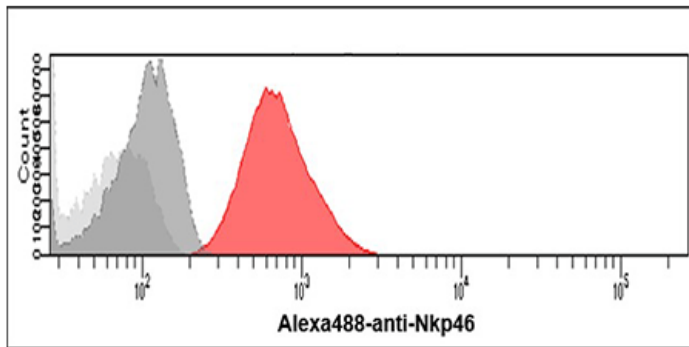
Ordering Information

[Order Online »](#)

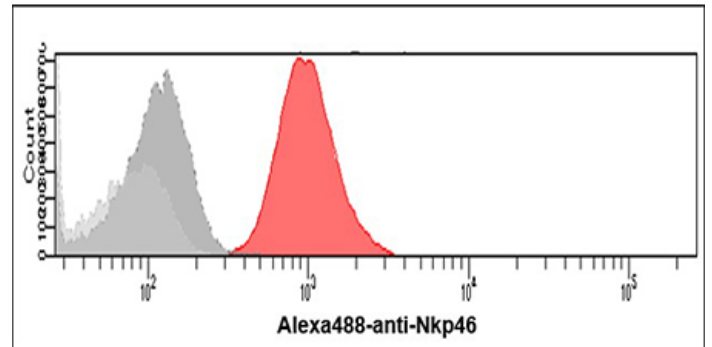
ENZ-ABS735-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of Nkp46/NCR1 in HeLa cells. The cell was stained at 2-5 μg for 1×10^6 cells (red). A Goat anti mouse IgG (Alexa Fluor 488) was used as the secondary antibody. Mouse monoclonal IgG was used as the isotype control (dark gray), cells without incubation with primary and secondary antibody was used as the negative control (light gray).



Flow cytometry analysis of Nkp46/NCR1 in PBMC. The cell was stained at 2-5 μg for 1×10^6 cells (red). A Goat anti mouse IgG (Alexa fluor 488) was used as the secondary antibody. Mouse monoclonal IgG was used as the isotype control (darkgray), cells without incubation with primary and secondary antibody was used as the negative control (light gray).

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Natural cytotoxicity triggering receptor 1 isoform a, NCR1, CD335
Application	ELISA, Flow Cytometry
Clone	n1D9
Formulation	Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.
Host	Mouse
Immunogen	Recombinant human Nkp46 (aa 22-255) purified from E. coli.
Isotype	IgG1κ
Purity Detail	Protein G affinity purified
Source	Purified from ascites.
Species Reactivity	Human
UniProt ID	O76036
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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