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 resion, 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-ABS736-0100 100μl

Manuals, SDS & CofA

View Online »

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/NCR1 (aa 22-255) purified

from E. coli.

Isotype IgG1ĸ

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human

UniProt ID 076036

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

Last modified: May 29, 2024



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

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

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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