

NKp30/NCR3 monoclonal antibody (AT38D9)

NCR3, also known as NKP30, is a natural cytotoxicity receptor (NCR) that may aid NK cells in the lysis of tumor cells. The encoded protein interacts with CD3-zeta (CD247), a T-cell receptor.

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

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 3, CD337

Application ELISA, WB

Clone AT38D9

Formulation Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.

Host Mouse

Immunogen Recombinant human NCR3 (aa 19-138) purified from E. coli.

Isotype IgG1κ

Purity Detail Protein A affinity purified

Source Purified from ascites.

Species Reactivity Human

UniProt ID O14931

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



ENZO LIFE SCIENCES,
Last modified **May 29, 2024**
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

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

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

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