

NKG2A polyclonal antibody

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

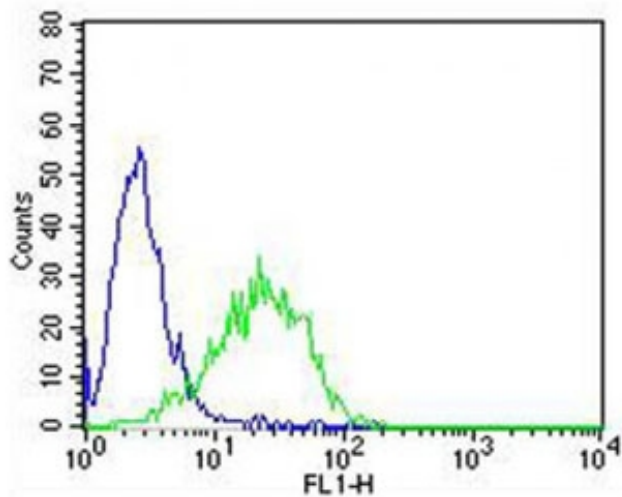
Ordering Information

[Order Online »](#)

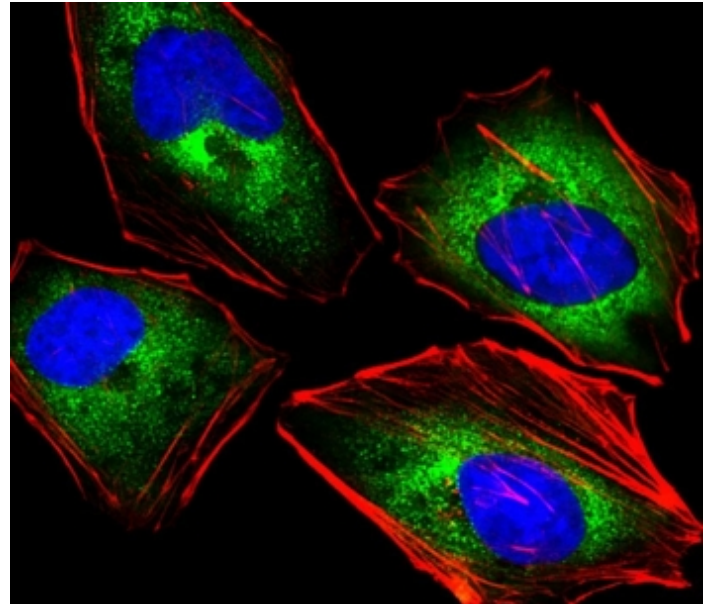
ENZ-ABS908-0200	0.2ml
-----------------	-------

Manuals, SDS & CofA

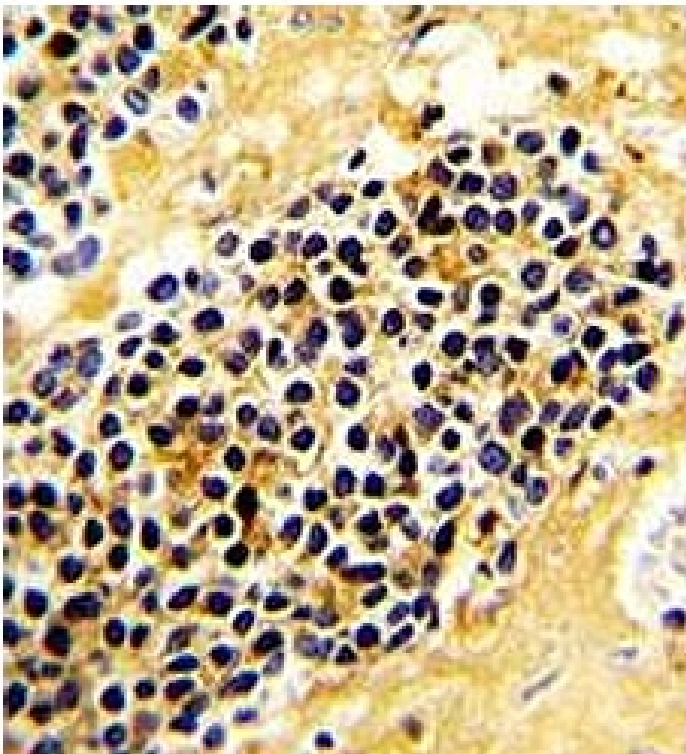
[View Online »](#)



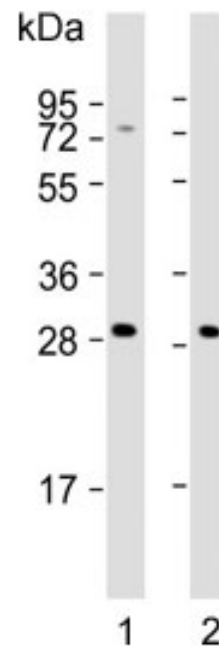
Flow cytometry testing of fixed and permeabilized human Jurkat cells with NKG2A antibody; Blue=isotype control, Green= NKG2A antibody.



Immunofluorescent staining of fixed and permeabilized human HeLa cells with NKG2A antibody (green), DAPI nuclear stain (blue) and anti-Actin (red).



IHC testing of FFPE human kidney carcinoma with NKG2A antibody. HIER: steam section in pH6 citrate buffer for 20 min and allow to cool prior to staining.



Western blot testing of human 1) K562 and 2) HepG2 cell lysate with NKG2A antibody. Predicted molecular weight: 26-40 kDa depending on glycosylation level.

Handling & Storage

Handling	Aliquot the NKG2A antibody and store frozen at -20°C or colder. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	KLRC1, NK cell receptor A, CD159a
Application	Flow Cytometry, IF, IHC (PS), WB
Formulation	Liquid. In 1X PBS, pH 7.4, containing 0.09% sodium azide.
Host	Rabbit
Immunogen	Human NKG2A (aa 180-206).
Isotype	Ig
Purity Detail	Antigen affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (1:25, 1×10 ⁶ cells)Immunofluorescence (1:25)Immunohistochemistry (FFPE) (1:25)Western blot (1:500-1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P26715
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)