## NKG2A polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

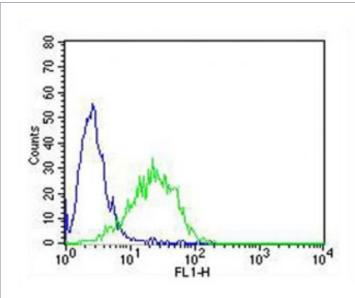
Order Online »

ENZ-ABS908-0200

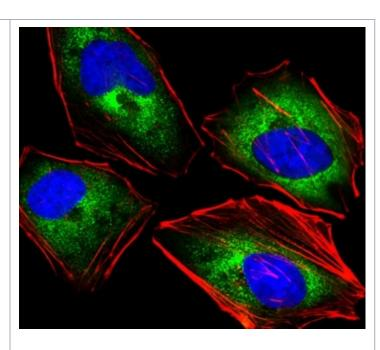
0.2ml

Manuals, SDS & CofA

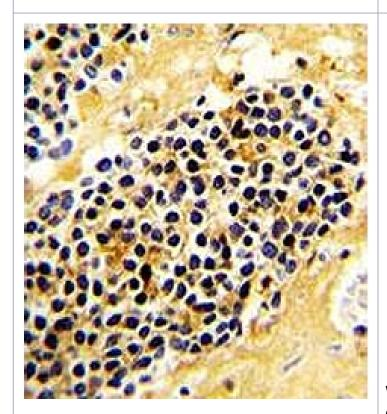
**View Online »** 



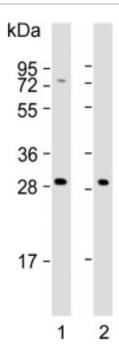
Flow cytometry testing of fixed and permeabilized human Jurkat cells with NKG2A antibody; Blue=isotype control, Green= NKG2A antibody.



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, 1×106 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

.



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