

# CD57 monoclonal antibody (NK-1)

Anti-CD57 marks a subset of lymphocytes known as natural killer (NK) cells. Follicular center cell lymphomas often contain many NK cells within the neoplastic follicles. Anti-CD57 also stains neuroendocrine cells and their derived tumors, including carcinoid tumor and medulloblastoma. Anti-CD57 can also be useful in separating type B3 thymoma from thymic carcinoma when combined with a panel that includes antibodies against GLUT1, CD5 and CEA.

CD57, also known as HNK-1 (human natural killer-1), is a cell surface marker associated with specific immune cell populations. It plays a crucial role in immune response modulation, neurobiology, and various pathological conditions. CD57 is a glycoprotein expressed on the surface of certain immune cells, including natural killer (NK) cells, T cells, and subsets of B cells. In the immune system, CD57 serves as a marker for identifying specific subsets of lymphocytes. Elevated CD57 expression on NK cells, for example, is linked to enhanced cytotoxicity and a more robust immune response against infected or aberrant cells. Additionally, CD57 can be used to distinguish memory T cell subsets, providing insights into the immune system's adaptive memory.

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

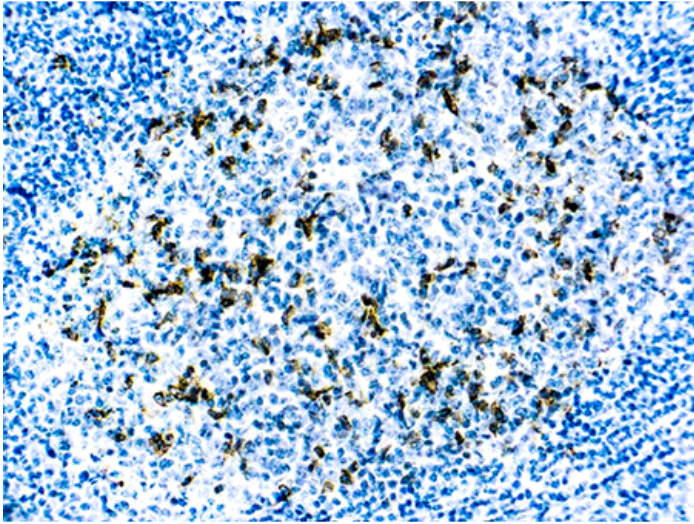
## Ordering Information

[Order Online »](#)

|                 |       |
|-----------------|-------|
| ENZ-ABS852-1000 | 1ml   |
| ENZ-ABS852-0100 | 100µl |

## Manuals, SDS & CofA

[View Online »](#)



Formalin fixed paraffin embedded human tonsil stained with CD57 antibody.

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## Product Details

Alternative Name Human natural killer-1, HNK-1

Application IHC (PS)

Application Notes This product is intended for qualitative immunohistochemistry with normal and neoplastic formalin-fixed, paraffin-embedded tissue sections, to be viewed by light microscopy.

Clone NK-1

Formulation This product is supplied as a purified immunoglobulin and contains less than 0.1% sodium azide as a preservative.

Host Mouse

Immunogen Human peripheral blood mononuclear cells.

Isotype IgM

Positive Control Spleen, Tonsil

Recommendation Dilutions/Conditions Antibody dilutions should be appropriately adjusted and verified according to the detection system used. Optimal conditions must be determined individually for each application.

Species Reactivity Human

### 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-  
be@enzolifesciences.com](mailto:info-be@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)

Last modified: May 29, 2024

For local distributors and detailed product information visit us online: [enzo.com](http://enzo.com) All product names, logos, and brands are the property of their respective owners.