

CD73 monoclonal antibody (AT89E9)

CD73 (Ecto-5-nucleotidase, 5-NT) is a membrane-bound glycoprotein that functions to hydrolyze extracellular nucleoside monophosphates enabling the uptake of adenosine, inosine, and guanosine into the cell and their subsequent reconversion into ATP and GTP. The expression of CD73 on neuronal cells has been described during developmental processes and, on lymphocytes, CD73 serves as a maturation marker, being absent from the surface of both immature B and T cells.

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

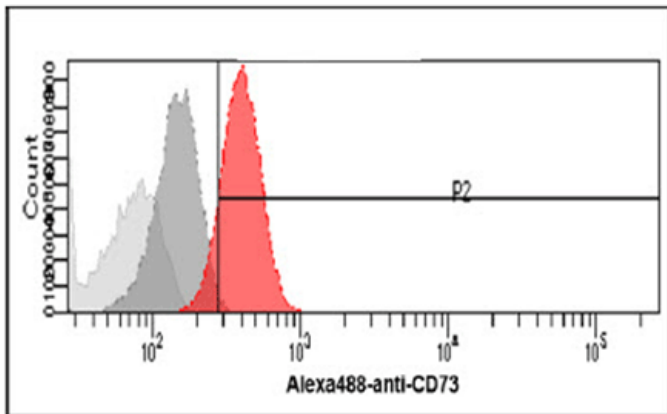
Ordering Information

[Order Online »](#)

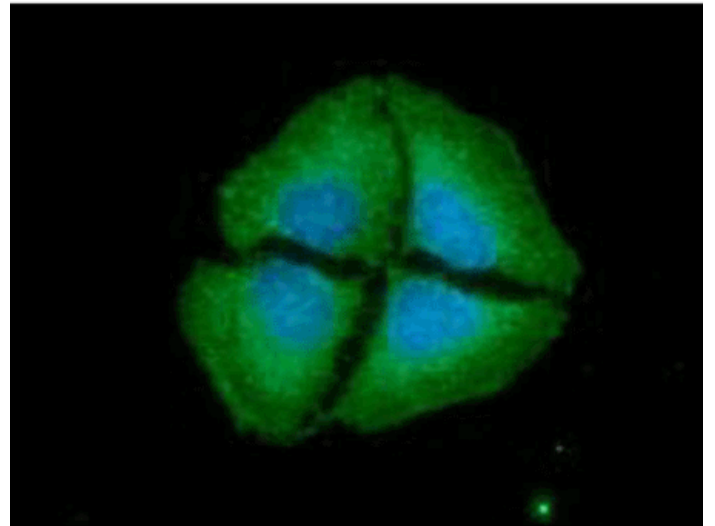
ENZ-ABS719-0100	100µl
-----------------	-------

Manuals, SDS & CofA

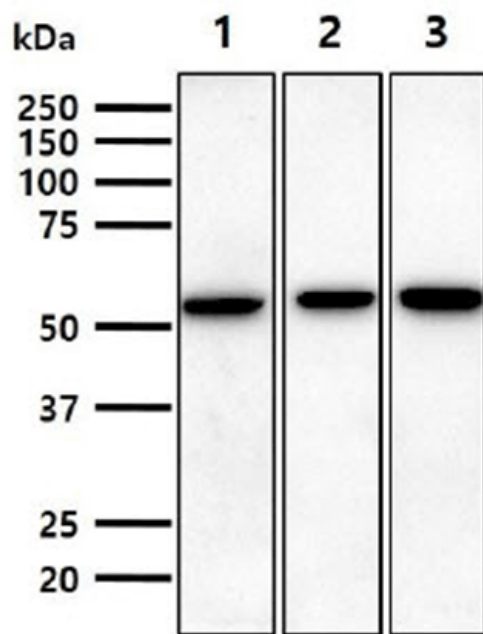
[View Online »](#)



Flow cytometry analysis of CD73 in A431 cells. The cell was stained with ENZ-ABS719 at 2-5 μ g for 1×10^6 cells (red). A Goat anti mouse IgG (Alexa fluor 488) was used as the secondary antibody. Mouse monoclonal IgG was used as the isotype control (dark gray), cells without incubation with primary and secondary antibody was used as the negative control (light gray).



ICC/IF analysis of CD73 in HeLa cells. The cell was stained with ENZ-ABS719 (1:100). The secondary antibody (green) was used Alexa Fluor 488. DAPI was stained the cell nucleus (blue).



The cell lysates (40 μ g) were resolved by SDS-PAGE, transferred to PVDF membrane and probed with anti-human CD73 antibody (1:1000). Proteins were visualized using a goat anti-mouse secondary antibody conjugated to HRP and an ECL detection system. Lane 1: HepG2 cell lysate. Lane 2: WiDr cell lysate. Lane 3: A431 cell lysate.

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 Ecto-5'-nucleotidase, 5'-NT

Application ELISA, Flow Cytometry, ICC, IF, WB

Clone AT89E9

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

Host Mouse

Immunogen Recombinant human CD73 (aa 27-252) purified from E. coli.

Isotype IgG2bk

Purity Detail Protein G affinity purified

Recommendation Dilutions/Conditions Western Blot (1:1,000). Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from ascites.

Species Reactivity Human

UniProt ID P21589

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)