

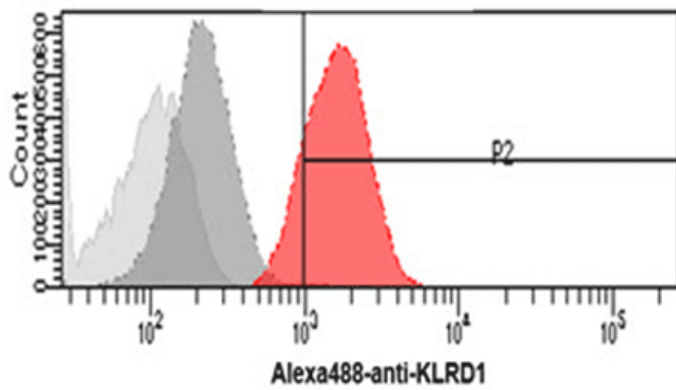
CD94 monoclonal antibody (AT13E3)

The protein encoded by CD94 gene is a lectin, cluster of differentiation and a receptor that is involved in cell signaling and is expressed on the surface of natural killer cells in the innate immune system. CD94 pairs with the NKG2 molecule as a heterodimer. The CD94/NKG2 complex, on the surface of natural killer cells interacts with Human Leukocyte Antigen (HLA)-E on target cells.

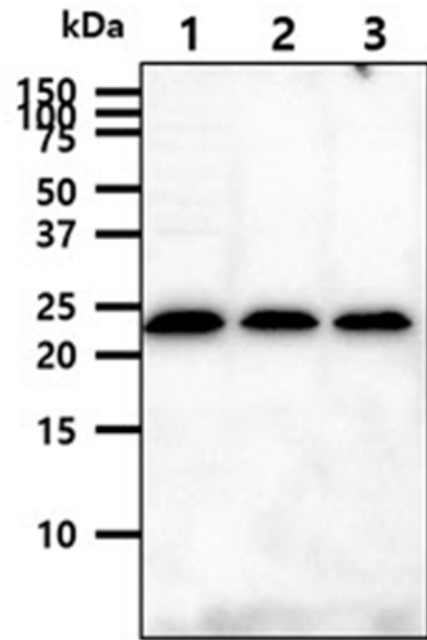
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of KLRD1 in PBMC cells. The cell was stained with ENZ-ABS723 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).



The cell lysates (40 μ g) were resolved by SDS-PAGE, transferred to PVDF membrane and probed with anti-human KLRD1 antibody (1:1000). Proteins were visualized using a goat anti-mouse secondary antibody conjugated to HRP and an ECL detection system. Lane 1: K562 cell lysate. Lane 2: MCF7 cell lysate. Lane 3: A549 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	Natural killer cells antigen CD94, KP43, Killer cell lectin-like receptor subfamily D member 1, NK cell receptor
Application	ELISA, Flow Cytometry, WB
Clone	AT13E3
Formulation	Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 10% glycerol.
Host	Mouse
Immunogen	Recombinant human KLRD1 (aa 32-179) purified from E. coli.
Isotype	IgG1k
Purity Detail	Protein A affinity purified
Source	Purified from ascites.
Species Reactivity	Human
UniProt ID	Q13241

Worry-free Guarantee

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



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

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

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

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

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

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