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.

Ordering Information

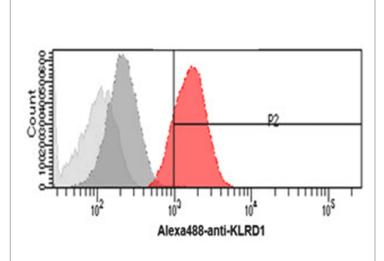
Order Online »

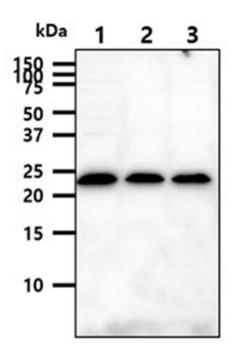
ENZ-ABS723-0100

100µl

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 1×10⁶ 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 (40µg) were resolved by SDS-PAGE, transferred to PVDF membrane and probed with antihuman 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 IgG1ĸ

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

