

# CD158i monoclonal antibody (5F2)

The KIR family consists of transmembrane glycoproteins of the Ig superfamily expressed on human Nk cells and a subset of human T cells which they are involved in recognition of either MHC class I molecules or unknown ligand on target cells and inhibit cytotoxic activities. KIR2DS4 is an activating receptor of KIR family.

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

## Ordering Information

[Order Online »](#)

ENZ-ABS730-0100	100µl
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## 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** Killer cell immunoglobulin-like receptor 2DS4, KIR2DS4, NKAT-8, Natural killer-associated transcript 8

**Application** ELISA, WB

**Clone** 5F2

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

**Host** Mouse

**Immunogen** Recombinant human KIR2DS4 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** P43632

**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)