

CD158b2 monoclonal antibody (190IIC311)

Reacts with extracellular domain of the KIR2DL3, a member of the Killer cell immunoglobulin-like receptor (KIR) family. 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.

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

Ordering Information

[Order Online »](#)

ENZ-ABS728-0100	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 2DL3

Application ELISA, ICC, IF, WB

Clone 190IIC311

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

Host Mouse

Immunogen Recombinant human KIR2DL3 (aa 19-161) purified from E. coli.

Isotype IgG2ak

Purity Detail Protein G affinity purified

Source Purified from ascites.

Species Reactivity Human

UniProt ID P43628

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)