

Anti-Human Interferon Lambda Receptor 1, Clone MMHLR-1 (MAb)

Binds to extracellular domains of human interferon lambda receptor.

Ordering Information

[Order Online »](#)

PBL-21885-1	50µg
-------------	------

Manuals, SDS & CofA

[View Online »](#)

- Neutralizes Human IFN Lambda Receptor 1
- Suitable for Flow Cytometry

Handling & Storage

Short Term Storage -20°C

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IFN Lambda Receptor 1

Application Flow Cytometry

Clone MMHLR-1

Formulation Phosphate – buffered saline (PBS) containing 0.1% bovine serum albumin (BSA)

Host Mouse

Immunogen Human IFN Lambda Receptor 1 extracellular domains

Isotype IgG1 kappa

Purity ≥ 95%

Species Reactivity Human

Specificity Binds to extracellular domains of human interferon lambda receptor



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-
be@enzolifesciences.com](mailto:info-be@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)