

Anti-Human Interleukin- 28A/Interferon Lambda 2, neutralizing, Clone

Neutralizes human IL-28A/interferon lambda 2

MMHL-2 (MA)

Ordering Information

[Order Online »](#)

PBL-21720-1

200µg

Manuals, SDS & CofA

[View Online »](#)

- Mouse monoclonal **antibody** against **Human IL-28A/IFNL2**
- **Neutralizes** human IL-28A/IFNL2
- ~3% and 8% cross-reactivity is observed with human IFNL1 and human IFNL3, respectively

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 IL-28A/IFNL2

Clone MMHL-2

Formulation Supplied frozen in phosphate-buffered saline (PBS) containing 0.1% bovine serum albumin (BSA).

Host Mouse

Isotype IgG1 kappa

Purity > 95%

Purity Detail Protein G affinity chromatography purified

Species Reactivity Human



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

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

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

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