

Anti-Human Interleukin- 28B/Interferon Lambda 3, neutralizing, Clone

Neutralizes human IL-28B/interferon lambda 3

MMHL-3 (MA)

Ordering Information

[Order Online »](#)

PBL-21730-1

100µg

Manuals, SDS & CofA

[View Online »](#)

- Mouse monoclonal **antibody** against Human IL-28B/IFNL3
- **Neutralizes** human IL-28B/IFNL3
- Less than 0.1% cross-reactivity is observed with human IFNL1 and human IFNL2

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-28B/IFNL3

Clone MMHL-3

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

Host Mouse

Immunogen Human IL28B/Interferon Lambda 3

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