

Anti-Mouse Interferon Beta, Clone RMMB-1 (MAb)

Binds to mouse interferon beta; weakly neutralizes mouse interferon beta

Ordering Information

[Order Online »](#)

PBL-22400-1	250µg
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	PBL-22400-1 (Unlabelled) should be shipped on dry ice. Storage -20°C. PBL-22400-3 (FITC-labelled) should be shipped on blue ice. Storage +4°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	IFN- β , Interferon- β
Clone	RMMB-1
Formulation	Liquid. In PBS containing 0.1% BSA.
Host	Rat
Immunogen	Mouse IFN- β (interferon- β).
Isotype	IgG1
Species Reactivity	Mouse



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