

Mouse Interleukin-28B/Interferon Lambda 3

IFN- λ 2, IFN- λ 3 and IFN- λ 1, also named IL-28A, IL-28B and IL-29, respectively, are newly identified class II cytokine receptor ligands that are distantly related to members of the IL-10 family and type I IFN family.

Ordering Information

[Order Online »](#)

PBL-12820-1	25 μ g
-------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	IL-28B/IFN-λ3
Endotoxin Content	< 1 EU/μg
Formulation	Lyophilized.
MW	~19.7kDa
Purity	≥95%
Source	Produced in <i>E. coli</i> .



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