

# Mouse Interleukin-28B/Interferon Lambda 3, carrier-free

IFN- $\lambda$ 2, IFN- $\lambda$ 3 and IFN- $\lambda$ 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-12821-1	25 $\mu$ 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](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)