

Human Interleukin-29/Interferon Lambda 1, carrier-free

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-11826-1 | 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-29/IFN-λ1

Formulation Lyophilized.

MW ~21.4kDa

Purity ≥95%

Source Produced in mouse myeloma cell line NSO.



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