

# Human Interferon Alpha B2 (Alpha 8)

## Ordering Information

[Order Online »](#)

PBL-11115-1	0.1MU
-------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Interferon- $\alpha$ B2, IFN- $\alpha$ 8, Interferon- $\alpha$ 8

**Endotoxin Content** < 1 EU/ $\mu$ g

**Formulation** Liquid. In PBS containing 0.1% BSA.

**MW** ~19.5kDa

**Purity**  $\geq$ 95%

**Purity Detail** Purified from *E. coli* by a combination of ion exchange, hydrophobic interaction and size exclusion chromatography

**Quantity** 0.1ml

**Source** Produced in *E. coli*.



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)