

# Human Interferon Alpha C (Alpha 10)

## Ordering Information

[Order Online »](#)

PBL-11120-1	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- $\alpha$ C, IFN- $\alpha$ 10, Interferon- $\alpha$ 10
Endotoxin Content	< 1 EU/ $\mu$ g
Formulation	Liquid. In PBS containing 0.1% BSA.
MW	~19.4kDa
Purity	$\geq$ 95%
Purity Detail	Purified from <i>E. coli</i> by a combination of ion exchange, hydrophobic interaction and size exclusion chromatography
Quantity	0.1ml
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)