

# Human Interferon Alpha I (Alpha 17)

## Ordering Information

[Order Online »](#)

|             |       |
|-------------|-------|
| PBL-11150-1 | 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- $\alpha$ 1, IFN- $\alpha$ 17, Interferon- $\alpha$ 17  |
| Endotoxin Content | < 1 EU/ $\mu$ g  |
| Formulation       | Liquid. In PBS containing 0.1% BSA.  |
| MW                | ~19.3kDa   |
| 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)