

**STD**

Human IFN alpha kit - Standard

Product reference :

**62HIFNACDA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 03A  
- Conditioning buffer\_\_\_\_\_ Phosphate 50 mM pH7.0 +  
BSA 1% + stabilizers  
- Appearance\_\_\_\_\_ Lyophilized form  
- Storage conditions\_\_\_\_\_ 2-8 °C  
- Expiration date\_\_\_\_\_ 2022.02.20

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Ab)	3 600.00 pg	N/A

**PLEASE NOTE**

Reconstitute with 0.300 mL H<sub>2</sub>O.

***Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)***

