

**IL2**

Human IL2 kit - Standard

Product reference :

**62HIL02CDA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 05A  
- Conditioning buffer\_\_\_\_\_ Phosphate 50 mM pH 7.0 +  
BSA 1% + stabilizers  
- Appearance\_\_\_\_\_ Lyophilized form  
- Storage conditions\_\_\_\_\_ 2-8 °C  
- Expiration date\_\_\_\_\_ 2022.12.12

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (MAb)	7 200.00 pg	15 000

*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)*

