

**STD**

Human IL1 beta kit - Standard

Product reference :

**62HIL1BCDA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 04A  
- Conditioning buffer\_\_\_\_\_ Phosphate 50 mM pH7.0 +  
BSA 1% + stabilizers  
- Appearance\_\_\_\_\_ Lyophilized form  
- Storage conditions\_\_\_\_\_ 2-8 °C  
- Expiration date\_\_\_\_\_ 2023.06.04

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Ab)	5 460.00 pg	N/A

*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)*