

## Batch information

### STD

Human IL1 beta kit - Standard

Product reference :

**62HIL1BCDA**

### Product characteristics :

- Lot Number	04B
- 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

### PLEASE NOTE

*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)*

