

## Batch information

**STD** Human IL23 kit - Standard

Product reference : **62HIL23CDA**

### 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	2024.06.02

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	27 000.00 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)*

