

## Batch information

**STD** HTRF Human IL17 kit - Std  
Product reference : **62HIL17CDA**

### Product characteristics :

- Lot Number	02A
- Conditioning buffer	Phosphate 50 mM pH7.0 + BSA 1% + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2025.07.18

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (IL)	10 500.00 pg	N/A

### PLEASE NOTE

Reconstitute with 0.350 mL H<sub>2</sub>O.

The expiration date of batch 02A has been extended to 2025-07-18 after validation by QC.

**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)**

