

Batch information

STD Human IL17 kit - Standard

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	2023.07.18

Weight indications :

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

PLEASE NOTE

Reconstitute with 0.350 mL H₂O.

The expiration date of batch 02A has been extended to 2023-07-18 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

