

Batch information

STD Human IL17 kit - Standard

Product reference : **62HIL17CDA**

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	2024.07.21

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (IL)	7 800.00 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

