

Batch information

IL5 HTRF Human IL5 kit - Std
Product reference : **62HIL05CDA**

Product characteristics :

- Lot Number	03A
- Conditioning buffer	Phosphate 50 mM pH 7.0 + BSA 1% + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2025.04.18

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (MAb)	7 500.00 pg	N/A

PLEASE NOTE

The expiration date of batch 03A has been extended to 2025-04-18 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

