

STD

Human IL1 alpha kit - Standard

Product reference :

62HIL1ACDA

Product characteristics :

- Lot Number _____ 02A
- Conditioning buffer _____ Phosphate 50 mM pH 7.0 +
BSA 1% + stabilizers
- Appearance _____ Lyophilized form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2023.03.28

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (MAb)	3 750.00 pg	15 000

PLEASE NOTE

The expiration date of batch 02A has been extended to 2023-03-28 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

