

STD

Human CCL5 kit - Standard

Product reference :

62HCCL5CDA

Product characteristics :

- Lot Number_____ 01A
- Conditioning buffer_____ Phosphate 50 mM pH7.0 +
BSA 1% + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2022.12.18

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (A)	1 800.00 pg	N/A

PLEASE NOTE

Reconstitute with 0.300 mL H₂O.

The expiration date of batch 01A has been extended to 2022-12- 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

