

STD

Human CCL2 kit - Standard

Product reference :

62HCCL2CDA**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_____ 2021.12.19

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (MAb)	2 250.00 pg	1

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com