

STD Human CCL4 kit - Standard
Product reference : **62HCCL4CDA**

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_____ 2022.03.13

Weight indications :

| | Quantity per vial | MW (Da) |
|---------------------|-------------------|---------|
| Active Moiety (MAb) | 4 500.00 pg | N/A |

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

