

LTC4 Standard
Product reference :

LTC4 kit calibrator
64LC4C0A

Product characteristics :

- Lot Number_____ 08A
- Conditioning buffer_____ ethanol
- Appearance_____ Frozen form
- Storage conditions_____ -60 °C
- Expiration date_____ 2022.03.27
- Volume per vial_____ 20 µL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (LC4)	2.00 µg	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