

EPIgeneous Lysis Buffer C

Product reference :

Epigeneous lysis buffer C - 500 tests

62EL3FDD

Product characteristics :

- Lot Number_____ 08A
- Conditioning buffer_____ Epigeneous lysis buffer C - 500 tests - Part 1 + Part 2
- Appearance_____ Frozen form
- Storage conditions_____ -20 °C
- Expiration date_____ 2024.01.08

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	500.00 tests	N/A

PLEASE NOTE

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