

**EPIgeneous Lysis Buffer C**

Product reference :

Epigeneous lysis buffer C - 10,000 tests

**62EL3FDH**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 09A  
- Conditioning buffer \_\_\_\_\_ Lysis buffer C - Part 1 + Part 2  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ -20 °C  
- Expiration date \_\_\_\_\_ 2025.03.17

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	10 000.00 tests	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](http://www.cisbio.com)*