

**Lysis buffer 2**

Product reference :

Lysis buffer 2 (4X) - 70 ml

**64LB2FDF**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 09A  
- Conditioning buffer \_\_\_\_\_ lysis buffer # 2  
- Appearance \_\_\_\_\_ Liquid form  
- Storage conditions \_\_\_\_\_ 2-8 °C  
- Expiration date \_\_\_\_\_ 2023.05.23  
- Volume per vial \_\_\_\_\_ 70 mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	70.00 mL	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)*