

**Lysis buffer # 4**  
Product reference :

Lysis buffer 4 - 130 ml  
**64LB4FDF**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 10A  
- Conditioning buffer \_\_\_\_\_ lysis buffer # 4  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ -20 °C  
- Expiration date \_\_\_\_\_ 2023.06.26  
- Volume per vial \_\_\_\_\_ 130 mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	130.00 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)*