

**lysis buf6**

Product reference :

Lysis & Detection Buffer 6 - 200 ml

**62CL6FDF**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 07A  
- Conditioning buffer\_\_\_\_\_ -  
- Appearance\_\_\_\_\_ Liquid form  
- Storage conditions\_\_\_\_\_ 2-8 °C  
- Expiration date\_\_\_\_\_ 2024.06.30

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	200.00 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)*