

**conj/lysis buf2**  
Product reference :

Lysis & Detection Buffer 2 - 200 ml  
**62CL2FDF**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 23A  
- Conditioning buffer \_\_\_\_\_ -  
- Appearance \_\_\_\_\_ Liquid form  
- Storage conditions \_\_\_\_\_ 2-8 °C  
- Expiration date \_\_\_\_\_ 2022.08.21

**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)*

