

**Buffer enz** Enzymatic buffer/kinase 5X - 10 ml  
Product reference : **62EZBFDC**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 12A  
- Conditioning buffer\_\_\_\_\_ -  
- Appearance\_\_\_\_\_ Liquid form  
- Storage conditions\_\_\_\_\_ 2-8 °C  
- Expiration date\_\_\_\_\_ 2023.12.13

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	10.00 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)*

