

**buffer 1**

Kinase Binding buffer - 200 mL

Product reference :

**62KBBRDF**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 01A  
- Conditioning buffer \_\_\_\_\_ -  
- Appearance \_\_\_\_\_ Liquid form  
- Storage conditions \_\_\_\_\_ 2-8 °C  
- Expiration date \_\_\_\_\_ 2024.01.17

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Kin)	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)*

