

**Cellular kinase lysis buffer#4**

Phospho-total protein lysis buffer #4 - 130 mL

Product reference :

**64KL4FDF**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 11A  
- Conditioning buffer\_\_\_\_\_ stock solution 4X  
- Appearance\_\_\_\_\_ Liquid form  
- Storage conditions\_\_\_\_\_ 2-8 °C  
- Expiration date\_\_\_\_\_ 2024.06.02

**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)*

