

**cell.phospho/total protein LB#3** Phospho-total protein lysis buffer #3 - 130 mL  
Product reference : **64KL3FDF**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 11A  
- Conditioning buffer \_\_\_\_\_ Stock solution 4X  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ 2-8 °C  
- Expiration date \_\_\_\_\_ 2024.01.10

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	130.00 mL	N/A

**PLEASE NOTE**

See corresponding package insert for use.

***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)***