

Cellular kinase lysis buffer#4

Phospho-total protein lysis buffer #4 - 130 mL

Product reference :

64KL4FDF

Product characteristics :

- Lot Number _____ 12A
- Conditioning buffer _____ stock solution 4X
- Appearance _____ Liquid form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2025.01.15

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

