

Cellular kinase lysis buffer

Product reference :

Phospho-total protein lysis buffer #1 - 130 mL

64KL1FDF**Product characteristics :**

- Lot Number	30A
- Conditioning buffer	Stock solution 4X
- Appearance	Liquid form
- Storage conditions	2-8 °C
- Expiration date	2023.11.27
- Volume per vial	130 mL

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