

Batch information

Cellular kinase lysis buffer

Phospho-total protein - Lysis buffer 1 (4X)- 130 mL

Product reference :

64KL1FDF**Product characteristics :**

- Lot Number	26A
- Conditioning buffer	Stock solution 4X
- Appearance	Liquid form
- Storage conditions	2-8 °C
- Expiration date	2022.07.27
- Volume per vial	130.00 mL

Weight indications :

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

PLEASE NOTE

The expiration date of batch 26A has been extended to 2022-07-27 after validation by QC.

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

