

## Batch information

### Cellular kinase lysis buffer

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

Product reference :

**64KL1FDF**

### Product characteristics :

- Lot Number	32A
- Conditioning buffer	Stock solution 4X
- Appearance	Liquid form
- Storage conditions	2-8 °C
- Expiration date	2026.02.18
- Volume per vial	130.00 mL

### Weight indications :

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

### PLEASE NOTE

*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)*

