

Cellular kinase lysis buffer

Product reference :

Phospho-total protein lysis buffer #1 - 130 mL

64KL1FDF

Product characteristics :

- Lot Number _____ 29A
- Conditioning buffer _____ Stock solution 4X
- Appearance _____ Liquid form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2022.12.12
- 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