

cell.phospho/total protein LB#3 Phospho-total protein lysis buffer #3 - 130 mL
Product reference : **64KL3FDF**

Product characteristics :

- Lot Number _____ 12A
- Conditioning buffer _____ Stock solution 4X
- Appearance _____ Frozen form
- Storage conditions _____ 2-8 °C
- Expiration date _____ 2024.02.21

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

