

Batch information

cell.phospho/total protein LB#3 HTRF P-T prot. - Lysis Buf.3 (4X) 130 mL

Product reference : **64KL3FDF**

Product characteristics :

- Lot Number	13A
- Conditioning buffer	Stock solution 4X
- Appearance	Liquid form
- Storage conditions	2-8 °C
- Expiration date	2025.12.21

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

