

Batch information

Cellular kinase lysis buffer

HTRF P-T prot. - Lysis Buf.1 (4X) 130 mL

Product reference :

64KL1FDF

Product characteristics :

- Lot Number	33A
- Conditioning buffer	Stock solution 4X
- Appearance	Liquid form
- Storage conditions	2-8 °C
- Expiration date	2027.02.02
- 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

