

Batch information

Cellular kinase lysis buffer#4 HTRF P-T prot. - Lysis Buf.4 (4X) 130 mL
Product reference : **64KL4FDF**

Product characteristics :

- Lot Number	13A
- Conditioning buffer	stock solution 4X
- Appearance	Liquid form
- Storage conditions	2-8° °C
- Expiration date	2026.10.11

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

