

## Batch information

**Cellular kinase lysis buffer**

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

Product reference :

**64KL1FDF****Product characteristics :**

- Lot Number	21A
- Conditioning buffer	Stock solution 4X
- Appearance	Liquid form
- Storage conditions	2-8 °C
- Expiration date	2022.08.09

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Buf)	130.00 mL	N/A

**PLEASE NOTE**

The expiration date of batch 21A has been extended to 2022-08-09 after validation by CQ.

***Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)***

