

Batch information

Cellular kinase blocking reagent Phospho-total protein - Blocking reagent- 6 mL

Product reference : **64KB1AAD**

Product characteristics :

- Lot Number	22A
- Conditioning buffer	Stock solution 100X
- Appearance	Liquid form
- Storage conditions	2-8 °C
- Expiration date	2023.07.03

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	6.00 mL	N/A

PLEASE NOTE

See corresponding package insert for use.

The expiration date of batch 22A has been extended to 2023-07-03 after validation by QC

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

