

Batch information

cell.phosp/tot protein BR

Phospho-total protein - Blocking reagent (100X)- 2 mL

Product reference :

64KB1AAC**Product characteristics :**

| | |
|-----------------------|---------------------|
| - Lot Number | 21B |
| - Conditioning buffer | Stock solution 100X |
| - Appearance | Frozen form |
| - Storage conditions | <-20 °C |
| - Expiration date | 2023.02.14 |

Weight indications :

| | Quantity per vial | MW (Da) |
|---------------------|-------------------|---------|
| Active Moiety (Buf) | 2.00 mL | N/A |

PLEASE NOTE

The expiration date of batch 21B has been extended to 2023-02-14 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

