

Batch information

cell.phosp/tot protein BR

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

Product reference :

64KB1AAC

Product characteristics :

- Lot Number	22C
- Conditioning buffer	Stock solution 100X
- Appearance	Frozen form
- Storage conditions	<-20 °C
- Expiration date	2023.07.03

Weight indications :

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

PLEASE NOTE

The expiration date of batch 22C 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

