

Batch information

cell.phosp/tot protein BR

Phospho-total protein - Blocking reagent- 2 ml

Product reference :

64KB1AAC

Product characteristics :

- Lot Number	25A
- Conditioning buffer	Stock solution 100X
- Appearance	Frozen form
- Storage conditions	<-20 °C
- Expiration date	2023.12.17

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Buf)	2.00 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

