

cell.phosp/tot protein BR

Phospho-total protein blocking reagent - 2 ml

Product reference :

64KB1AAC

Product characteristics :

- Lot Number _____ 23F
- Conditioning buffer _____ Stock solution 100X
- Appearance _____ Frozen form
- Storage conditions _____ <-20 °C
- Expiration date _____ 2022.09.19

Weight indications :

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

PLEASE NOTE

See corresponding package insert for use.

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

