

Batch information

cell.phosp/tot protein BR

HTRF P-T prot. - Block. reag.(100X) 2 mL

Product reference :

64KB1AAC**Product characteristics :**

- Lot Number	23F
- Conditioning buffer	Stock solution 100X
- Appearance	Frozen form
- Storage conditions	<-20 °C
- Expiration date	2024.09.19

Weight indications :

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

PLEASE NOTE

The expiration date of batch 23F has been extended to 2024-09-19 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

