

Batch information

Cellular kinase blocking reagent HTRF P-T prot. - Block. reag.(100X) 2 mL

Product reference : **64KB1AAC**

Product characteristics :

- Lot Number	27A
- Conditioning buffer	Stock solution 100X
- Appearance	Frozen form
- Storage conditions	<-20 °C
- Expiration date	2023.06.20

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

