

## Batch information

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

Product reference : **64KB1AAC**

### Product characteristics :

- Lot Number	23C
- Conditioning buffer	Stock solution 100X
- Appearance	Frozen form
- Storage conditions	<-20 °C
- Expiration date	2023.09.19

### Weight indications :

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

### PLEASE NOTE

The expiration date of batch 23C has been extended to 2023-09-19 after validation byQC

***Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)***

