

## Batch information

**cell.phosp/tot protein BR**

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

Product reference :

**64KB1AAC**

### Product characteristics :

- Lot Number	26A
- Conditioning buffer	Stock solution 100X
- Appearance	Frozen form
- Storage conditions	<-20 °C
- Expiration date	2025.02.24

### 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](http://www.cisbio.com)*

