

**Tracer** Staurosporine-Red - 1 nmol  
**Product reference :** **62KB01REDC**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 02A  
- Conditioning buffer \_\_\_\_\_ -  
- Appearance \_\_\_\_\_ Frozen form  
- Storage conditions \_\_\_\_\_ <-16 °C  
- Expiration date \_\_\_\_\_ 2023.02.19

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Kin)	40.00 µl	N/A

**PLEASE NOTE**

The expiration date of batch 02A has been extended to 2023-02-19 after validation

***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)***

