

**Tracer**

Staurosporine-Red - 20 nmol

Product reference :

**62KB01REDE**

**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) | 800.00 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)***

