

**Control lysate**

STING total kit control lysate

Product reference :

**64NTGTDA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 03A  
- Conditioning buffer\_\_\_\_\_ Phospho total protein lysis buffer  
- Appearance\_\_\_\_\_ Frozen form  
- Storage conditions\_\_\_\_\_ -60 °C  
- Expiration date\_\_\_\_\_ 2022.12.10  
- Volume per vial\_\_\_\_\_ 150 µL  
- Active Moiety Concentration (Ab)\_\_\_\_\_ 1 X

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Ab)	150.00 µl	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)***

