

**Control lysate**

CREB phospho-S133 kit control lysate

Product reference :

**64CRETDA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 13A  
- Conditioning buffer\_\_\_\_\_ Cellular kinase lysis buffer  
- Appearance\_\_\_\_\_ Frozen form  
- Storage conditions\_\_\_\_\_ -60 °C  
- Expiration date\_\_\_\_\_ 2022.06.19  
- Volume per vial\_\_\_\_\_ 150 µL  
- Active Moiety Concentration (CL)\_\_\_\_\_ 1 X

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (CL)	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)***

