

**Control lysate**

GSK3 alpha phospho-S21 kit control lysate

Product reference :

**64GPATDA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 001  
- Conditioning buffer\_\_\_\_\_ Lysis Buffer  
- Appearance\_\_\_\_\_ Frozen form  
- Storage conditions\_\_\_\_\_ -60 °C  
- Expiration date\_\_\_\_\_ 2022.09.19  
- Volume per vial\_\_\_\_\_ 150 µL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Ab)	150.00 µl	N/A

**PLEASE NOTE**

The expiration date of batch 01 has been extended to 2022-09-19 after validation by QC.

***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)***