

**Control lysate**

Alpha-synuclein phospho-S129 control lysate

Product reference :

**6FSYNTDA**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 08A  
- Conditioning buffer\_\_\_\_\_ Lysis Buffer  
- Appearance\_\_\_\_\_ Frozen form  
- Storage conditions\_\_\_\_\_ -60 °C  
- Expiration date\_\_\_\_\_ 2022.07.17  
- Volume per vial\_\_\_\_\_ 150 µL

**Weight indications :**

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

*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)*