

Batch information

Control lysate Alpha-synuclein total kit - Standard
Product reference : **6FNSYCDA**

Product characteristics :

- Lot Number 06A
- Conditioning buffer Lysis Buffer
- Appearance Lyophilized form
- Storage conditions 2-8 °C
- Expiration date 2024.03.12
- Volume per vial 300.00 µL
- Active Moiety Concentration (Ab) 500.00 ng/mL

Weight indications :

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

PLEASE NOTE

The expiration date of batch 06A has been extended to 2024-03-12 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

