

Batch information

Control lysate

HTRF GSK3 alpha p-S21 kit - Ctrl lysate

Product reference :

64GPATDA

Product characteristics :

- Lot Number		1
- Conditioning buffer	Lysis Buffer	
- Appearance	Frozen form	
- Storage conditions	-60 °C	
- Expiration date	2024.09.19	
- Volume per vial	150.00 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 2024-09-19 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

