

Control lysate

GSK3 beta phospho-S9 kit control lysate

Product reference :

64GPBTDA**Product characteristics :**

- Lot Number	04A
- Conditioning buffer	Cellular kinase lysis buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2022.08.09
- 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 04A has been extended to 2022-08-09 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

