

Batch information

Control lysate HTRF GSK3 beta p-S9 kit - Ctrl lysate
Product reference : **64GPBTDA**

Product characteristics :

- Lot Number	03A
- Conditioning buffer	Cellular kinase lysis buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2023.08.18
- Volume per vial	150.00 µL
- Active Moiety Concentration (Ab)	1.00 X

Weight indications :

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

PLEASE NOTE

The expiration date of batch 03A has been extended to 2023-08-18 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

