

## Batch information

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

### Product characteristics :

- Lot Number	04A
- Conditioning buffer	Cellular kinase lysis buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.08.09
- 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 04A has been extended to 2024-08-09 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](http://www.cisbio.com)**

