

## Batch information

### Control lysate

GAPDH Housekeeping kit - control lysate

Product reference :

**64GAPDHTDA**

### Product characteristics :

- Lot Number	02A
- Conditioning buffer	Lysis Buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.08.10
- Volume per vial	150.00 µL
- Active Moiety Concentration (CL)	1.00 X

### Weight indications :

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

### PLEASE NOTE

*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)*

