

## Batch information

### Control lysate

HTRF b-Catenin total kit - Ctrl lysate

Product reference :

**64NCATDA**

### Product characteristics :

- Lot Number	02A
- Conditioning buffer	Lysis Buffer
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2024.05.17
- 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 02A has been extended to 2024-05-17 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)***

