

## Batch information

### control lysate

HTRF B-arrestin 2 total kit - Ctrl lys.

Product reference :

**64BAR2TTDA**

### Product characteristics :

- Lot Number	01A
- Conditioning buffer	lysis buffer #4 sans BR
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2023.11.25

### Weight indications :

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

*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)*

