

## Batch information

**Total H3 Cellular Control Lysate**      HTRF Total H3 cell. assay - Ctrl lysate  
Product reference :                      **62NH3TDA**

### Product characteristics :

- Lot Number	13A
- Conditioning buffer	Mix Lysis Buffer A
- Appearance	Frozen form
- Storage conditions	-60 °C
- Expiration date	2025.06.01
- Volume per vial	150.00 µL

### Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (CTL)	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)*

