

Total H3 Cellular Control Lysate Histone H3 total kit control lysate
Product reference : **64NH3TDA**

Product characteristics :

- Lot Number_____ 07A
- Conditioning buffer_____ Mix Lysis Buffer A
- Appearance_____ Frozen form
- Storage conditions_____ -60 °C
- Expiration date_____ 2022.01.21
- Volume per vial_____ 150 µL
- Active Moiety Concentration (CTL)_____ 1 X

Weight indications :

| | Quantity per vial | MW (Da) |
|---------------------|-------------------|---------|
| Active Moiety (CTL) | 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