

Control lysate

HIF-1 alpha kit control lysate

Product reference :

64HIFTDA

Product characteristics :

- Lot Number_____ 02A
- Conditioning buffer_____ lysis buffer
- Appearance_____ Frozen form
- Storage conditions_____ -60 °C
- Expiration date_____ 2022.08.08
- Volume per vial_____ 150 µL

Weight indications :

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

