

anti 6HIS-d2

Product reference :

MAb Anti 6HIS-d2 - 4 mg - CUSTOM LABELING

61HISDAJ**Product characteristics :**

| | |
|-------------------------------------|---------------------------------------|
| - Lot Number | 23B |
| - Packaging Unit | 200 000 tests/vial |
| - Conditioning buffer | PO4 0.1M pH7 + 0.1% BSA + stabilizers |
| - Appearance | Frozen form |
| - Storage conditions | <-16 °C |
| - Expiration date | 2029.05.18 |
| - Volume per vial | 8 mL |
| - Active Moiety Concentration (HIS) | 500 µg/mL |

Weight indications :

| | Quantity per vial | MW (Da) |
|---------------------|-------------------|---------|
| Active Moiety (HIS) | 4.00 mg | 150 000 |

Recommended quantity per well* :

| | in 384 well low volume plate (20µL) |
|----------------------|-------------------------------------|
| Active Moiety (HIS) | 20.00 ng |
| Number of tests/vial | 200 000 |

*Actual amount per well should be optimized according to assay conditions

PLEASE NOTE

The performance of assays using 6HIS-tag can be enhanced by incubating those at 4-8°C.

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com