

**anti 6HIS-d2**

MAb Anti 6HIS-d2 - 20,000 tests

Product reference :

**61HISDLB****Product characteristics :**

- Lot Number	24A
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	PO4 0.1M pH7 + 0.1% BSA + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2023.01.08
- Reconstitute each vial with	1 mL distilled H2O
- Active Moiety Concentration (HIS)	400 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (HIS)	400.00 µg	150 000

**Recommended quantity per well\* :**

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

\*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](http://www.cisbio.com)**