

anti 6HIS-d2
Product reference :

MAb Anti 6HIS-d2 - 4 mg - CUSTOM LABELING
61HISDAJ

Product characteristics :

- Lot Number_____ 23C
 - 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_____ 2029.07.05
 - 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