

MAB6HIS-XL
Product reference :

MAB Anti 6HIS-XL665 - 5,000 tests
61HISXLA

Product characteristics :

- Lot Number_____ 62A
- Packaging Unit_____ 5 000 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1% BSA + Stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2022.09.26
- Reconstitute each vial with_____ 0.25 mL distilled H2O
- Active Moiety Concentration (MAB)_____ 400 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (MAB)	100.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (MAB)	20.00 ng
Number of tests/vial	5 000

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

PLEASE NOTE

After reconstitution the concentration of Active Moiety (MAB) is 400 µg/mL.

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