

anti 6HIS-d2
Product reference :

MAb Anti 6HIS-d2 - 4 mg
61HISDAJ

Product characteristics :

- Lot Number_____ 19A
- Conditioning buffer_____ PO4 0.1M pH7 + 0.1% BSA
- Appearance_____ Frozen form
- Storage conditions_____ <-16 °C
- Expiration date_____ 2026.06.07
- 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

PLEASE NOTE

The performance of assays using 6HIS-tag can be enhanced by incubating those at 2-8°C.

Further information as well as a specific application note regarding conjugate-related calculations are available on our HTRF web site: www.htrf.com/products