

**anti 6HIS-d2**

Product reference :

**61HISDAJ**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 16A
- Packaging Unit \_\_\_\_\_ 200 000 tests/vial
- Conditioning buffer \_\_\_\_\_ PO4 0.1M pH7 + 0.1% BSA
- Appearance \_\_\_\_\_ Frozen form
- Storage conditions \_\_\_\_\_ <-16 °C
- Expiration date \_\_\_\_\_ 2024.12.04
- 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 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](http://www.htrf.com/products)**