

Batch information

anti 6HIS-d2

HTRF MAb Anti 6HIS-d2 - 1K pts

Product reference :

61HISDLF

Product characteristics :

- Lot Number	23D
- Packaging Unit	1 000 tests/vial
- Conditioning buffer	PO4 0.1M pH7 + 0.1% BSA + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2022.12.02
- Reconstitute each vial with	0.25 mL distilled H2O
- Active Moiety Concentration (HIS)	80.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (HIS)	20.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (HIS)	20.00 ng
Number of tests/vial	1 000

*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

