

Batch information

MAB 6HIS-Tb

HTRF MAb Anti 6HIS-Tb - 1K pts

Product reference :

61HISTLF

Product characteristics :

- Lot Number	13A
- Packaging Unit	1 000 tests/vial
- Conditioning buffer	PO4 100 mM pH7 + stabilizers + BSA 0.5%
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2025.01.05
- Reconstitute each vial with	0.25 mL distilled H2O
- Active Moiety Concentration (MAb)	8.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (MAb)	2.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (MAb)	2.00 ng
Number of tests/vial	1 000

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

PLEASE NOTE

The use of KF is not compulsory in HTRF assays using Lumi4-Tb conjugate.

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

