

Batch information

anti-GST-Tb

MAb Anti GST-Tb cryptate - 20,000 tests

Product reference :

61GSTTLB

Product characteristics :

- Lot Number	10A
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.5% BSA+ Stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8°C °C
- Expiration date	2023.09.19
- Reconstitute each vial with	1 mL distilled H2O
- Active Moiety Concentration (Mab)	50.00 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Mab)	50.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	2.50 ng
Number of tests/vial	20 000

*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 expiration date of batch 10A has been extended to 2023-09-19 after validation by QC.

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com

