

anti-GST-Tb

MAb Anti GST-Tb cryptate - 1,000 tests

Product reference :

61GSTTLF

Product characteristics :

- Lot Number_____ 09A
- Packaging Unit_____ 1 000 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.5% BSA+
Stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2023.01.04
- Reconstitute each vial with_____ 0.25 mL distilled H2O
- Active Moiety Concentration (Mab)_____ 10 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Mab)	2.50 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	2.50 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 expiration date of batch 09A has been extended to 2023-01-04 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