

anti-GST-Tb

MAb Anti GST-Tb cryptate - 1 mg

Product reference :

61GSTTAH

Product characteristics :

- Lot Number_____	10B
- Packaging Unit_____	400 000 tests/vial
- Conditioning buffer_____	Phosphate 0.1M pH7 + 0.5% BSA+ stabilizers
- Appearance_____	Frozen form
- Storage conditions_____	<-16 °C
- Expiration date_____	2029.02.11
- Volume per vial_____	4 mL
- Active Moiety Concentration (Mab)_____	250 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Mab)	1.00 mg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	2.50 ng
Number of tests/vial	400 000

*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.

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