

anti-GST-Tb

MAb Anti GST-Tb cryptate - 1 mg

Product reference :

61GSTTAH

Product characteristics :

- Lot Number_____ 11A
- 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_____ 2029.05.12
- 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