

Mab anti MBP-Tb
Product reference :MAB Anti MBP-Tb cryptate - 5,000 tests
61MBPTAA**Product characteristics :**

- Lot Number_____ 02A
- Packaging Unit_____ 5 000 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1% BSA +
0.1% Tween
- Appearance_____ Frozen form
- Storage conditions_____ <-16 °C
- Expiration date_____ 2027.05.09
- Volume per vial_____ 0.25 mL
- Active Moiety Concentration (Ab)_____ 30 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Ab)	7.50 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	1.50 ng
Number of tests/vial	5 000

*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