

**Mab anti MBP-Tb**  
Product reference :

MAb Anti MBP-Tb cryptate - 20,000 tests  
**61MBPTAB**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 02A  
- Packaging Unit\_\_\_\_\_ 20 000 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\_\_\_\_\_ 1 mL  
- Active Moiety Concentration (Ab)\_\_\_\_\_ 30 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Ab)	30.00 µg	150 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (Ab)	1.50 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.

**Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: [www.cisbio.com](http://www.cisbio.com)**