

Batch information

Mab anti MBP-Tb

MAB Anti MBP-Tb cryptate - 5,000 tests

Product reference :

61MBPTAA

Product characteristics :

- Lot Number	02B
- 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)	30.00 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

