

**MABFLAG M2-Tb**  
Product reference :MAB Anti FLAG M2-Tb cryptate - 1 mg - CUSTOM LABELING  
**61FG2TAH****Product characteristics :**

- Lot Number	18A
- Packaging Unit	769 230 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1% de BSA + 0.1% de Tween
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2029.05.18
- 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)	1.30 ng
Number of tests/vial	769 230

\*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)***