

MABFLAG M2-Tb
Product reference :

MAb Anti FLAG M2-Tb cryptate - 1 mg - CUSTOM LABELING
61FG2TAH

Product characteristics :

- Lot Number	14A
- Packaging Unit	769 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.1% de BSA + 0.1% de Tween
- Appearance	Frozen form
- Storage conditions	-20 °C
- Expiration date	2025.07.16
- Volume per vial	1.575 mL
- Active Moiety Concentration (Mab)	635 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	769 000

*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 HTRF web site: www.htrf.com/products