

MABFLAG M2-XL
Product reference :MAB Anti FLAG M2-XL665 - 20,000 tests
61FG2XLB**Product characteristics :**

- Lot Number_____ 26A
- Packaging Unit_____ 20 000 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1% BSA + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2022.09.19
- Reconstitute each vial with_____ 1 mL distilled H2O
- Active Moiety Concentration (Mab)_____ 0.4 mg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Mab)	400.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	20.00 ng
Number of tests/vial	20 000

*Actual amount per well should be optimized according to assay conditions

PLEASE NOTE

The expiration date of batch 26A has been extended to 2022-09-19 after validation by QC

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com