

Anti FlagM2-d2
Product reference :MAb Anti FLAG M2-d2 - 1,000 tests
61FG2DLF**Product characteristics :**

- Lot Number_____ 12B
- Packaging Unit_____ 1 000 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1% BSA+ stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2022.08.29
- Reconstitute each vial with_____ 0.25 mL distilled H2O
- Active Moiety Concentration (Mab)_____ 80 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (Mab)	20.00 µg	150 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	20.00 ng
Number of tests/vial	1 000

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

PLEASE NOTE

The expiration date of batch 12A has been extended to 2022-08-29 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