

**Anti FlagM2-d2**  
Product reference :MAb anti Flag M2-d2 - 20,000 tests  
**61FG2DLB****Product characteristics :**

- Lot Number\_\_\_\_\_ 15A  
- 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\_\_\_\_\_ 2023.08.20  
- Reconstitute each vial with\_\_\_\_\_ 1 mL distilled H2O  
- Active Moiety Concentration (Mab)\_\_\_\_\_ 400 µg/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

**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)**