

**Anti FlagM2-d2**  
Product reference :

Monoclonal Anti-Flag M2 Antibody labeled with d2  
**61FG2DAJ**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 06A
- Packaging Unit \_\_\_\_\_ 200 000 tests/vial
- Conditioning buffer \_\_\_\_\_ Phosphate 0.1M pH7 + 0.1% BSA
- Appearance \_\_\_\_\_ Frozen form
- Storage conditions \_\_\_\_\_ -20 °C
- Expiration date \_\_\_\_\_ 2022.07.03
- Volume per vial \_\_\_\_\_ 8 mL
- Active Moiety Concentration (Mab) \_\_\_\_\_ 500 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (Mab)	4.00 mg	150 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (Mab)	20.00 ng
Number of tests/vial	200 000

\*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 HTRF web site: [www.htrf.com/products](http://www.htrf.com/products)**