



## Technical Data Sheet

**Anti FlagM2-d2**

HTRF MAb Anti FLAG-d2 - 20K pts

Product reference :

**61FGBDLB****Product characteristics :**

- Lot Number	01AR
- 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	2026.12.03
- Reconstitute each vial with	1 mL distilled H2O
- Active Moiety Concentration (Mab)	400.00 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

**PLEASE NOTE**

*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)*