

**61KAPDAB**

MAb Anti Human Kappa-d2 -20,000 tests - CUSTOM LABELING

Product reference :

**61KAPDAB**

**Product characteristics :**

- Lot Number \_\_\_\_\_ 007  
 - Packaging Unit \_\_\_\_\_ 59 938 tests/vial  
 - Conditioning buffer \_\_\_\_\_ Phosphate 0.1M pH7 + 0.1 % BSA  
 - Appearance \_\_\_\_\_ Frozen form  
 - Storage conditions \_\_\_\_\_ -80 °C  
 - Expiration date \_\_\_\_\_ 2028.10.07  
 - Volume per vial \_\_\_\_\_ 1.8 mL  
 - Active Moiety Concentration (Ab) \_\_\_\_\_ 666 µg/mL

**Weight indications :**

|                    | Quantity per vial | MW (Da)   |
|--------------------|-------------------|-----------|
| Active Moiety (Ab) | 1 198 756.00      | ng 50 000 |

**Recommended quantity per well\* :**

|                      | in 384 well<br>low volume<br>plate (20µL) |
|----------------------|---|
| Active Moiety (Ab)   | 20.00 ng                                  |
| Number of tests/vial | 59 938                                    |

\*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)**