

**MAb MBP-D2**  
Product reference :

Mouse monoclonal Ab anti Maltose binding protein labeled with d2  
**61MBPDAB**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 05A
- Packaging Unit\_\_\_\_\_ 20 000 tests/vial
- Conditioning buffer\_\_\_\_\_ Phosphate 0.1M pH7 + 0.1% BSA
- Appearance\_\_\_\_\_ Frozen form
- Storage conditions\_\_\_\_\_ <-16 °C
- Expiration date\_\_\_\_\_ 2023.10.06
- Volume per vial\_\_\_\_\_ 1.6 mL
- Active Moiety Concentration (Mab)\_\_\_\_\_ 250 µg/mL

**Weight indications :**

|                     | Quantity per vial | MW (Da) |
|---------------------|-------------------|---------|
| Active Moiety (Mab) | 0.40 mg           | 150 000 |

**Recommended quantity per well\* :**

|                      | in 384 well<br>low volume<br>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 HTRF web site: [www.htrf.com/products](http://www.htrf.com/products)**