

Batch information

SA-d2 Streptavidin-d2 - 20,000 tests
Product reference : **610SADLB**

Product characteristics :

| | |
|------------------------------------|--|
| - Lot Number | 23A |
| - 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 | 2024.08.03 |
| - Reconstitute each vial with | 1 mL distilled H2O |
| - Active Moiety Concentration (SA) | 1000.00 µg/mL |

Weight indications :

| | Quantity per vial | MW (Da) |
|--------------------|-------------------|---------|
| Active Moiety (SA) | 1 000.00 µg | 60 000 |

Recommended quantity per well* :

| | in 384 well low volume plate (20µL) |
|----------------------|-------------------------------------|
| Active Moiety (SA) | 50.00 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

