

**SA-d2**

Streptavidin-d2 - 20,000 tests

Product reference :

**610SADLB**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 20A
- 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\_\_\_\_\_ 2022.01.30
- Reconstitute each vial with\_\_\_\_\_ 1 mL distilled H2O
- Active Moiety Concentration (SA)\_\_\_\_\_ 1000 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (SA)	1.00 mg	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

**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)**