

SA-d2

Streptavidin-d2 - 1,000 tests

Product reference :

610SADLF

Product characteristics :

- Lot Number_____ 21A
- Packaging Unit_____ 1 000 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7+ 0.1% BSA + stabilizers
- Appearance_____ Lyophilized form
- Storage conditions_____ 2-8 °C
- Expiration date_____ 2023.06.22
- Reconstitute each vial with_____ 0.25 mL distilled H2O
- Active Moiety Concentration (SA)_____ 200 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (SA)	50.00 µg	60 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (SA)	50.00 ng
Number of tests/vial	1 000

*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