

## Batch information

**SA-d2** Streptavidin-d2 - 1,000 tests  
Product reference : **610SADLF**

### Product characteristics :

- Lot Number	23A
- 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	2024.08.03
- Reconstitute each vial with	0.25 mL distilled H2O
- Active Moiety Concentration (SA)	200.00 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

### 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](http://www.cisbio.com)**

