

**SA-d2**

Streptavidin-d2 - 5 mg

Product reference :

**610SADAC****Product characteristics :**

- Lot Number	19F
- Packaging Unit	100 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7+ 0.1% BSA + stabilizers
- Appearance	Frozen form
- Storage conditions	<-16 °C
- Expiration date	2028.11.26
- Volume per vial	5 mL
- Active Moiety Concentration (SA)	1 mg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (SA)	5.00 mg	60 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (SA)	50.00 ng
Number of tests/vial	100 000

\*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)**