

SA-d2

Streptavidin-d2 - 5 mg

Product reference :

610SADAC

Product characteristics :

- Lot Number_____ 15A
- Packaging Unit_____ 100 000 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7+ 0.1% BSA + stabilizers
- Appearance_____ Frozen form
- Storage conditions_____ -20 °C
- Expiration date_____ 2025.04.13
- Volume per vial_____ 5 mL
- Active Moiety Concentration (SA)_____ 1000 µg/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 HTRF web site: www.htrf.com/products