

SA-Tb

Streptavidin-Tb cryptate - 20,000 tests

Product reference :

610SATLB**Product characteristics :**

- Lot Number	08B
- Packaging Unit	20 000 tests/vial
- Conditioning buffer	Phosphate 0.1M pH7 + 0.5 % BSA + stabilizers
- Appearance	Lyophilized form
- Storage conditions	2-8 °C
- Expiration date	2022.12.18
- Reconstitute each vial with	1 mL distilled H2O
- Active Moiety Concentration (SA)	12 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (SA)	12.00 µg	60 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (SA)	0.60 ng
Number of tests/vial	20 000

*Actual amount per well should be optimized according to assay conditions

PLEASE NOTE

The use of KF is not compulsory in HTRF assays using Lumi4-Tb conjugate.

Further information as well as a specific application note regarding conjugate-related calculations are available on our Life Sciences web site: www.cisbio.com