

**SA-Tb**

Streptavidin-Tb cryptate - 1,000 tests

Product reference :

**610SATLF**

**Product characteristics :**

- Lot Number\_\_\_\_\_ 08B  
 - Packaging Unit\_\_\_\_\_ 1 000 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\_\_\_\_\_ 0.25 mL distilled H2O  
 - Active Moiety Concentration (SA)\_\_\_\_\_ 2.4 µg/mL

**Weight indications :**

	Quantity per vial	MW (Da)
Active Moiety (SA)	0.60 µg	60 000

**Recommended quantity per well\* :**

	in 384 well low volume plate (20µL)
Active Moiety (SA)	0.60 ng
Number of tests/vial	1 000

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