

SA-Tb

Streptavidin-Tb cryptate - 1 mg - CUSTOM LABELING
610SATAH

Product reference :

Product characteristics :

- Lot Number_____ 07A
- Packaging Unit_____ 1 666 666 tests/vial
- Conditioning buffer_____ Phosphate 0.1M pH7 + 0.1 % BSA + 0.1% tween
- Appearance_____ Frozen form
- Storage conditions_____ -20 °C
- Expiration date_____ 2025.04.14
- Volume per vial_____ 1.391 mL
- Active Moiety Concentration (SA)_____ 719 µg/mL

Weight indications :

	Quantity per vial	MW (Da)
Active Moiety (SA)	1.00 mg	60 000

Recommended quantity per well* :

	in 384 well low volume plate (20µL)
Active Moiety (SA)	0.60 ng
Number of tests/vial	1 666 666

*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 HTRF web site: www.htrf.com/products