

HTRF® Research Reagents

## Streptavidin-Tb Quality Control

Product number : 610SATLF Lot Number : 21RB

Manufacturing date : 07/12/2022 Document version : 1

### EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing Anti GST-D2 concentrations.

SA-Tb is lyophilized in PO4 buffer 100 mM pH7 ; 0.5% BSA ; Stabilizers.

All the reagents involved were diluted in PPI-Terbium detection buffer (#61DB10RDF).

### ASSAY FORMAT

SA-Tb + Pept-GST-biot. + Anti GST-D2

### REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
Anti GST-d2	15A	0.1 - 0.375 - 0.75 - 1.5 nMf
SA-Tb reference	14A	0.333 nMf
Pept-GST-Biot	CQ09	2.5 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
SA-Tb current batch	21RB	0.333 nMf

*Other reagents were the same as those used for the reference test.*

### EXPERIMENTAL PROCEDURE

Reconstitute SA-Tb vial with 250 µL H2O (if packaged under 5,000 tests or 1,000 tests) or 1 mL H2O (if packaged under 20,000 tests).

Reagents were dispensed as follows :

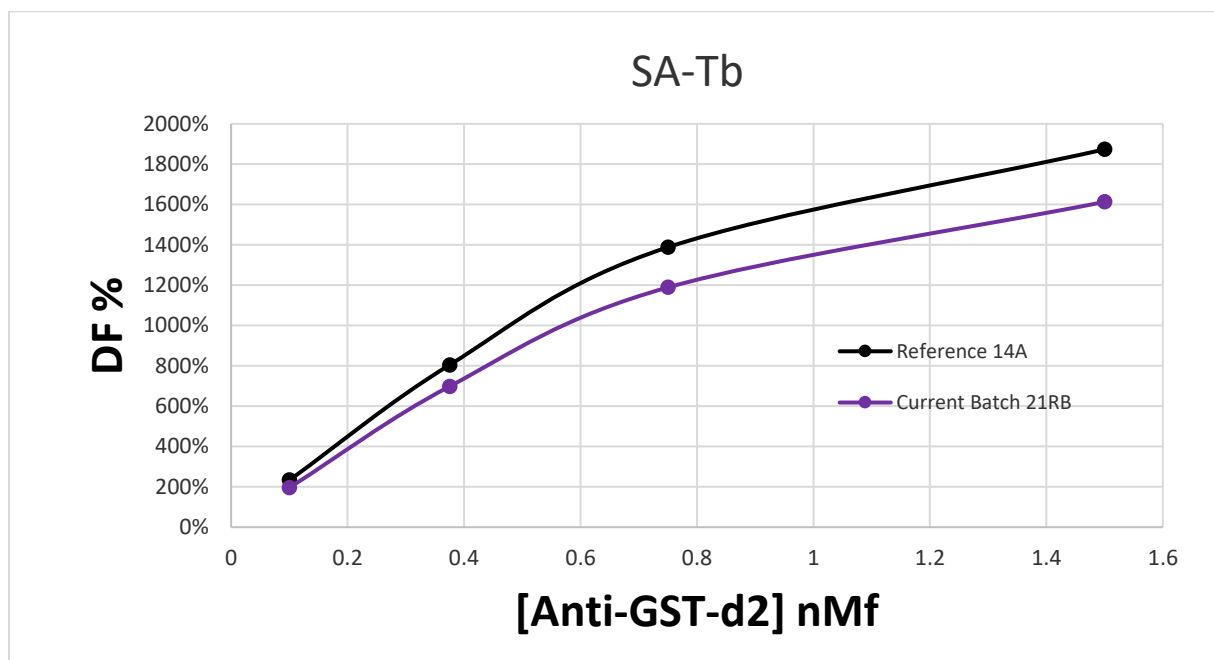
- 5 µL Pept-GST-Biotin (5 µL diluent for negative control)
- 10 µL D2 conjugate
- 5 µL Tb conjugate

Incubation took place for 120 min at room temperature. The plate was read on Phera Star ® under standard conditions.

## RESULTS

[Anti GST-D2] nMf		SA-Tb Reference	SA-Tb CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
0.1	% Delta F % CV	233 % 0.6 %	196% 1.4 %	-16.1%	-14.4%
0.375	% Delta F % CV	804 % 0.3%	698 % 0.3 %	-13.3%	
0.75	% Delta F % CV	1388 % 1.4 %	1189% 1.2%	-14.4%	± 20 %
1.5	% Delta F % CV	1874 % 1.0 %	1613 % 1.5%	-13.9%	

For the batch release, variation in delta F must be less than 20 %.



SA-Tb batch # 21RB

Accepted



Rejected



Operator's initials: MT

Date : 23/12/2022

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

revvity