

HTRF® Research Reagents

Anti-Mouse-Tb Quality Control

Product number : 61PAMTAF Lot Number : 03RC

Manufacturing date : 17/10/2022 Document version : 1

EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing conjugate XL concentrations (IgG1-XL).

Pab Anti Mouse IgG-Tb is frozen in Hepes buffer 50 mM pH 7 ; 0.1% BSA.

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

ASSAY FORMAT

Anti-Mouse-IgG-Tb or Anti-Rabbit IgG-Tb + IgG1-XL

REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
IgG1- XL	CQ01	0.21 - 3.33 nMf
Anti-Mouse-Tb	03B	62.5 ng/mLf
Anti-Rabbit-Tb	01A	75 ng/mLf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti-Mouse-Tb current batch	03RC	62.5 ng/mLf

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

EXPERIMENTAL PROCEDURE

Reagents were dispensed as follows :

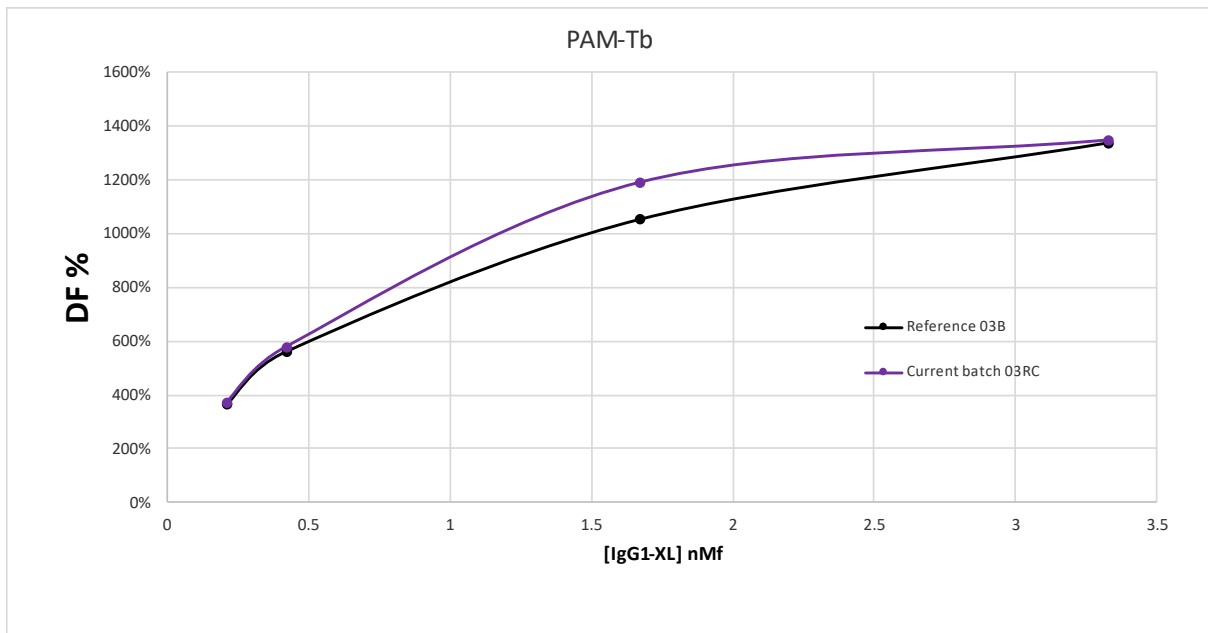
- 5 µL Tb conjugate
- 10 µL PPI - Terbium detection buffer
- 5 µL IgG1-XL

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

RESULTS

[IgG1-XL] nM _f		Anti-Mouse-Tb Reference	Anti-Mouse-Tb CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
0.21	% Delta F % CV	364 % 7.4 %	371% 0.5 %	2.11 %	4.71 %
0.42	% Delta F % CV	561 % 2.9 %	578 % 8.6 %	2.89 %	
1.67	% Delta F % CV	1 053 % 12.5 %	1 190 % 0.2 %	13.05 %	± 20 %
3.33	% Delta F % CV	1 338 % 4.6 %	1 348 % 2.1 %	0.81 %	

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



Anti-Mouse-Tb batch # 03RC

Accepted

Rejected

Operator's initials: ML

Date: 17/10/2022

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.

