

HTRF® Research Reagents

Mab anti MBP-Tb Quality Control

Product number : 61MPBTAA Lot Number : 03RB

Manufacturing date : 12/05/2023 Document version : 1

EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing Peptide concentrations (MBP-Biotin).

Mab anti MBP-Tb is frozen in buffer PO4 100 mM pH7; 0.1% BSA.

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

ASSAY FORMAT

SA-XL + MBP-Biotin + anti MBP-Tb

REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
SA-XL	178 A	0.37 – 1.1 – 3.3 – 10 nMf
Anti MBP-Tb _{Ref}	03A (61MBPTAB)	0.5 nMf
MBP-Biotin	02	0.37 – 1.1 – 3.3 - 10 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti MBP-Tb	03RB	0.5 nMf

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

EXPERIMENTAL PROCEDURE

Reagents were dispensed as follows :

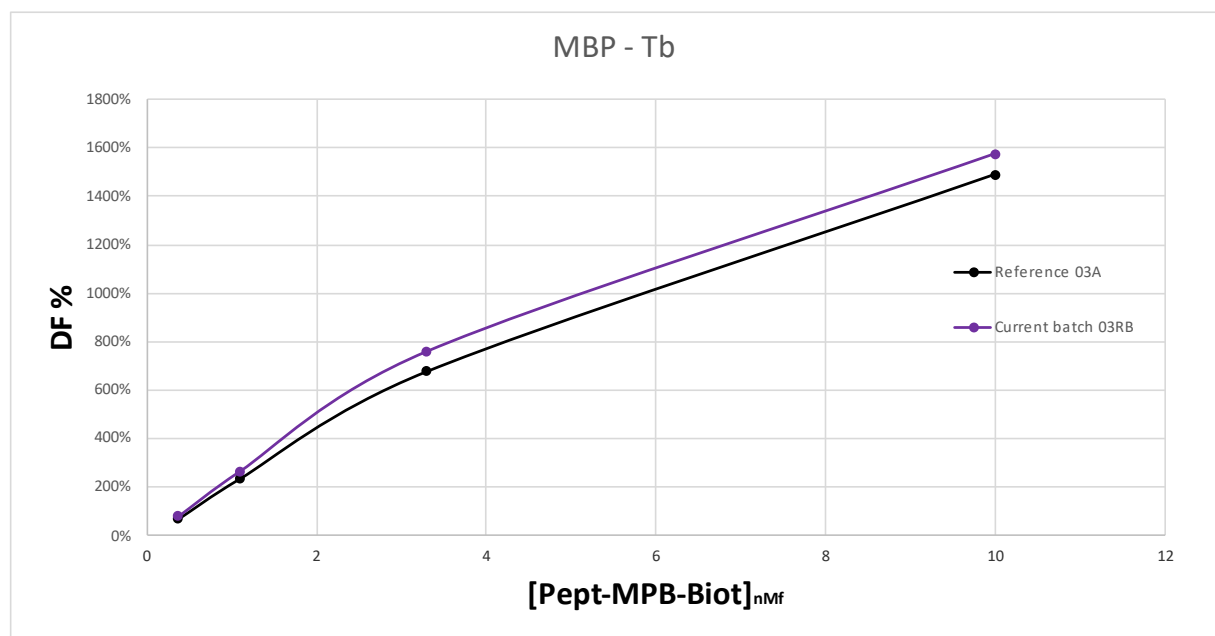
- 5 µl MBP-Biotin
- 10 µl XL conjugate
- 5 µl Tb Conjugate

Incubation took place for 180 min at room temperature. The plate was read on Phera Star ® under standard conditions.

RESULTS

[MBP Biotin] nMf		Anti MBP-Tb Ref Lot 03A	Anti MBP-Tb CQ 03RB	Variation	Mean Acceptable variation
0.37	% Delta F % CV	67 % 1.3%	77 % 1.7 %	15.1%	± 20 %
1.1	% Delta F % CV	234% 1.3%	264 % 2.4 %	12.8%	11.4%
3.3	% Delta F % CV	676 % 1.4 %	758 % 2.7%	12.2%	
10	% Delta F % CV	1488 % 1.5%	1574 % 1.9 %	5.7%	

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



MBP-Tb batch # 03RB

Accepted

Rejected

Operator's initials: ML

Date : 12/05/2023

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.

