

HTRF® Research Reagents

Anti MBP-K Quality Control

Product number : 61MBPKAA

Lot Number : 02RB

Manufacturing date : 30/01/2023

Document version : 1

EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing MBP-biotin concentrations.

Anti MBP-K is frozen in PO4 buffer 100 mM pH7 ; 0.1% BSA.

All the reagents involved were diluted in PPI-Europium detection buffer (#61DB9RDF).

ASSAY FORMAT

SA-XL + MBP-Biotin + Anti MBP-K

REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
SA-XL	170A	0.37 – 1.1 – 3.3 – 10 nMf
Anti MBP-K Reference	02A	0.67 nMf
MBP-Biotin	02	0.37 – 1.1 – 3.3 – 10 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti MBP-K current batch	02RB	0.67 nMf

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

EXPERIMENTAL PROCEDURE

Reagents were dispensed as follows :

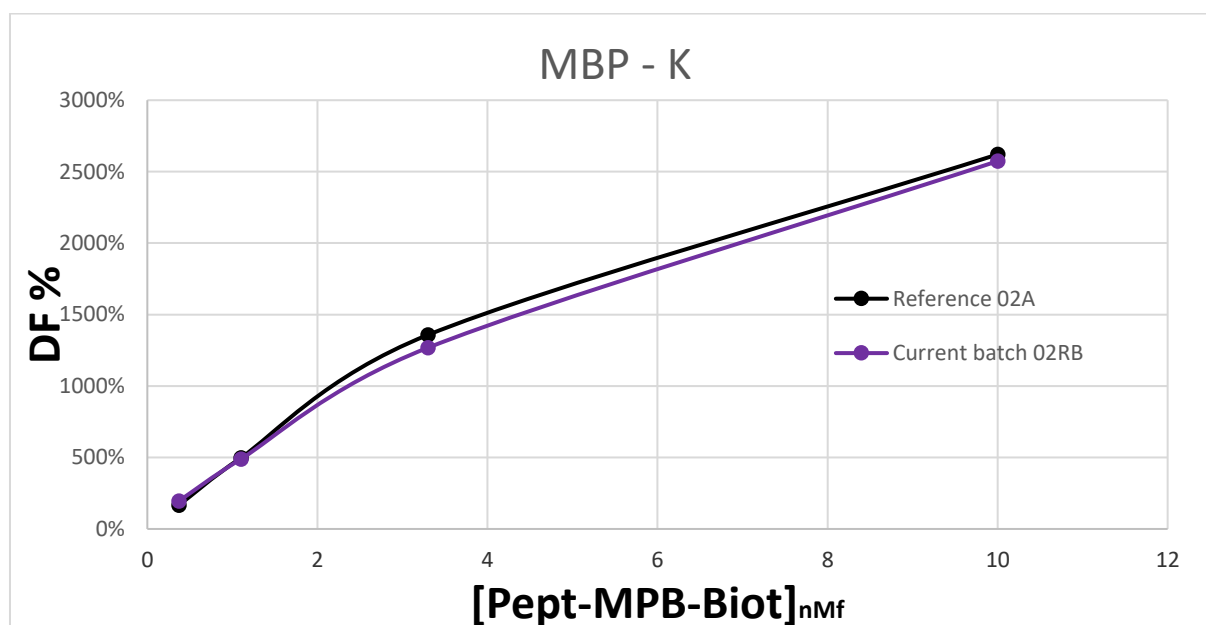
- 5 µL MBP-Biotin (5 µL diluent for negative control)
- 10 µL XL conjugate
- 5 µL K conjugate

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-K Reference	Anti MBP-K CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
0.37	% Delta F % CV	166 % 1.5%	195% 1.6 %	17.9 %	2.0%
1.1	% Delta F % CV	497 % -	489 % 2.6 %	-1.7 %	
3.3	% Delta F % CV	1358 % 4.4%	1268 % 3.2 %	-6.6 %	± 20 %
10	% Delta F % CV	2620% 4.5%	2573 % 0.2 %	-1.8%	

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



Anti MBP-K batch # 02RB

Accepted



Rejected



Operator's initials: MT

Date: 31/01/2023

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.

revvity