

HTRF® Research Reagents

Anti MBP-d2 Quality Control

Product number : 61MBPDAA Lot Number : 06CR

Manufacturing date : 26/10/2023 Document version : 1

EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing antibody concentrations (anti-MBP-d2).

Anti-MBP-d2 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

Anti-MBP-d2 + MBP-Biotin + Streptavidin-Cryptate

REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
Streptavidin-K	CQ02	2.33 nMf
Anti-MBP-d2 Ref	06A	0.1 - 0.5 - 1 - 2 - 4 nMf
MBP-Biotin	002	4.66 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti-MBP-d2 current batch	06CR	0.1 - 0.5 - 1 - 2 - 4 nMf

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

If applicable, streptavidin / biotin ratio was set up at : 1/2

EXPERIMENTAL PROCEDURE

Reagents were dispensed as follows :

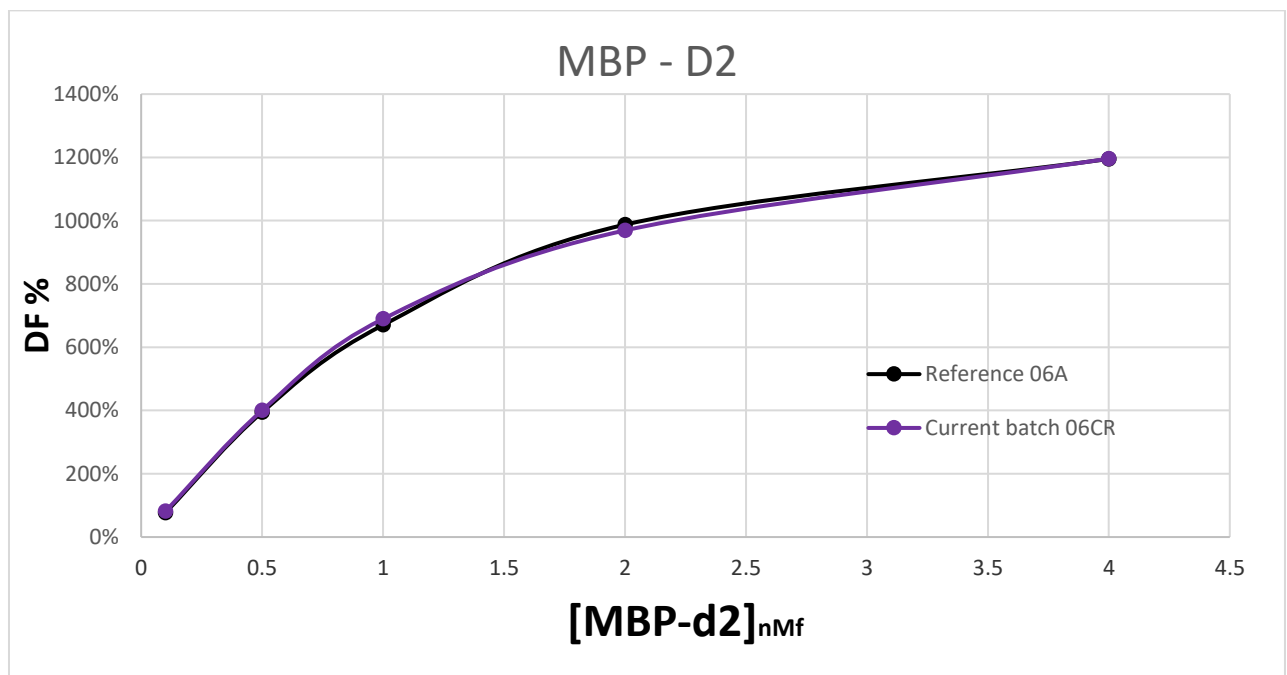
- 5 µL MBP-Biotin (5 µL diluent for negative control)
- 10 µL D2 conjugate
- 5 µL K conjugate

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

RESULTS

[Anti-MBP-d2] nM _f		Anti-MBP-d2 Ref	Anti-MBP-d2 CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
0.1	% Delta F % CV	77 % 1.3 %	81 % 1.6 %	6.0 %	1.7 %
0.5	% Delta F % CV	395 % 1.9 %	400 % 1.0 %	1.4 %	
1	% Delta F % CV	670 % 1.5 %	689 % 2.8 %	2.9 %	± 20 %
2	% Delta F % CV	987 % 4.8 %	969 % 0.9 %	-1.9 %	
4	% Delta F % CV	1195 % 1.4 %	1196 % 0.5 %	0.1 %	

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



Anti MBP-d2 batch # 06CR

Accepted



Rejected



Operator's initials: AL

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