

HTRF® Research Reagents

## Anti phospho p53-K Quality Control

Product number : 61P08KAZ Lot Number : 10RA

Manufacturing date : 07/01/2022 Document version : 1

### EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing Pept-PO4-biotin (n° 864) concentrations and at decreasing Pept-nonPO4-biotin concentrations (n° 876).

Anti phospho p53-K is frozen in HEPES buffer 50 mM pH 7, 0.1% BSA.

All the Xlent! conjugate and K conjugate involved were diluted in PPI Europium buffer #61DB9RDF

Peptide Biotin-PO<sub>4</sub> (n° 864) and Peptide Biotin- non PO<sub>4</sub> (n° 876) involved were diluted in PPI Europium buffer #61DB9RDF

### ASSAY FORMAT

Anti phospho p53-K + Peptide Biotin-PO<sub>4</sub> and Peptide Biotin- non PO<sub>4</sub> + SA-XL<sup>ent!</sup>

### REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
SA-XL <sup>ent!</sup>	25B	5 nMf
Anti phospho p53-k	09A	0.833 nMf
Pept-PO4-Biotin (n° 864)	CQ01	10 nMf (gradient de 0% à 100%)
Pept-PO4- non Biotin (n° 876)	CQ01	10 nMf (gradient de 0% à 100%)

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti phospho p53-k	10RA	0.833 nMf

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

### EXPERIMENTAL PROCEDURE

Reagents were dispensed as follows :

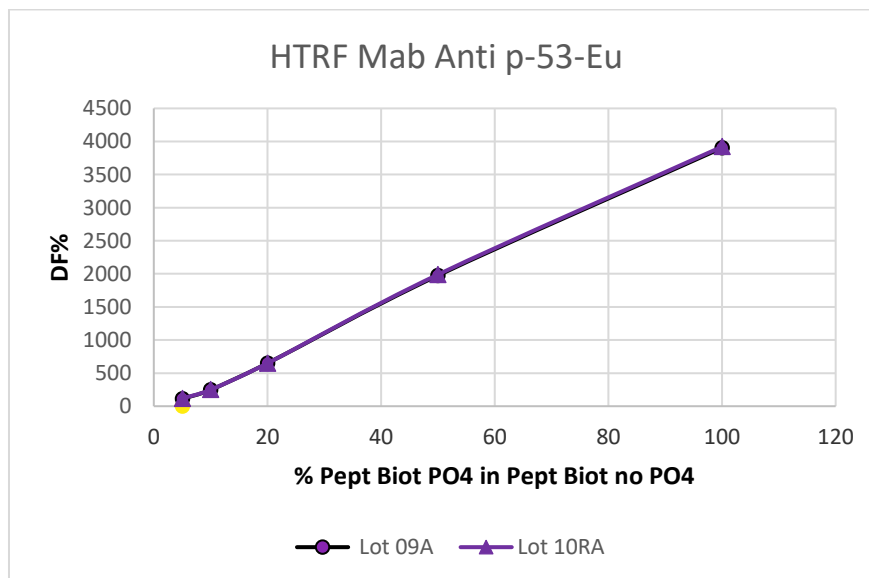
- 10 µL Pept-PO4-Biotin and Pept-non PO4-Biotin
- 5 µL XL conjugate
- 5 µL K conjugate

Incubation took place for 150 min at room temperature. The plate was read on Phera Star <sup>®</sup> under standard conditions.

## RESULTS

[Pept-PO4-Biotin] %		Anti phospho p53-k Ref	Anti phospho p53-k CURRENT BATCH	VARIATION	ACCEPTABLE VARIATION
5	% Delta F % CV	111 % 1.6 %	117 % 0.15 %	5.4 %	1.4 %
10	% Delta F % CV	248 % 1 %	250 % 1.4 %	0.8 %	
20	% Delta F % CV	649 % 1.1 %	649 % 2.5 %	0%	
50	% Delta F % CV	1975 % 0.2 %	1989 % 0.7 %	0.7%	± 20 %
100	% Delta F % CV	3903 % 2.1 %	3921 % 1.3 %	0.4 %	

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



Anti phospho p53-K batch #  
10RA

Accepted

Rejected

Operator's initials: ML

Date :13/01/2022

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