

Anti phospho p53-K Quality Control

Product reference : 61P08KAE

Batch # : 08C

1- Experimental conditions

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

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 50 mM Hepes pH 7, 0.4 M KF + 0.1 % BSA.

Peptide Biotin-PO₄ (n° 864) and Peptide Biotin- non PO₄ (n° 876) involved were diluted in Tp Hepes 50 mM pH=7 + 0.4 M KF + 0.1 % BSA.

2- Assay format

Anti phospho p53-K + Peptide Biotin-PO₄ and Peptide Biotin- non PO₄ + SA-XL^{ent!}

3- Reagents

	Reference lot number	Assay Concentration
SA-XL ^{ent!}	24A	5 nM _f
Anti phospho p53-K	08A	0.833 nM _f
Peptide Biotin-PO ₄ (n° 864)	CQ01	10 nM _f (gradient de 0% à 100%)
Peptide Biotin- non PO ₄ (n° 876)	CQ01	10 nM _f (gradient de 100% à 0%)
	Current batch lot number	Assay Concentration
Anti phospho p53-K	08C	0.833 nM _f

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

If applicable, Biotin / Streptavidin ratio was set up at : 2

4- Experimental procedure

Reagents were dispensed as follows :

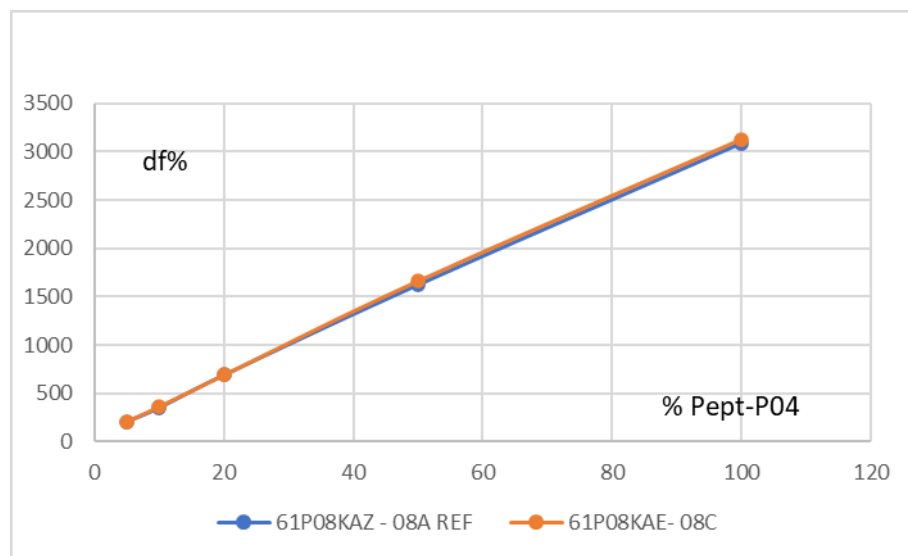
- 50 µl Peptide Biotin-PO₄ and Peptide Biotin- non PO₄
- 25 µl XL^{ent!} conjugate
- 25 µl K conjugate

Incubation took place for 150 min at room temperature. The plate was read on PHERA Star under standard conditions.

5- Results

Peptide Biot-PO ₄ (n°864) %		Anti phospho p53-K-08A Réf	Anti phospho p53-K-08C CQ	Variation / Réf	Variation Moyenne acceptable
5	% Delta F % CV	205 % 2.32%	208 % 0.71 %	1.86%	1.74 % ± 20 % CV < 10 %
10	% Delta F % CV	350 % 2.92%	362% 1.53 %	3.39%	
20	% Delta F % CV	688% 2.04 %	688% 1.41%	-0.10 %	
50	% Delta F % CV	1622 % 5.3 %	1660% 3.47%	2.35%	
100	% Delta F % CV	3089% 6.3%	3126% 0.02 %	1.17%	

The batch release , variation in delta F must be less than 20%



**Anti phospho p53-K batch # 08C
accepted**

Operator's initials : MB

Supervisor's initials and date
16/06/2020 LN