

HTRF® Research Reagents

## Anti phospho IκBa-K Quality Control

Product number : 61P06KAE Lot Number : 03RB

Manufacturing date : 04/08/2005 Document version : 1

### EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing Peptide Biotin-PO4 (n° 839) concentrations and at decreasing Peptide Biotin- non PO4 (n° 833) concentrations.

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

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

### ASSAY FORMAT

Anti phospho IκBa -K + Peptide Biotin-PO4 and Peptide Biotin-non PO4 + SA-XL<sup>ent!</sup>

### REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
SA-XL ent!	04F	5 nMf
Anti phospho IκBa -K reference	02F	31 ng/mLf
Peptide Biotin-PO4 (n° 839)	CQ01	10 nMf (gradient de 0% à 100%)
Peptide Biotin-non PO4 (n° 833)	CQ01	10 nMf (gradient de 100% à 0%)

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti phospho IκBa -K current batch	03RB	34 ng/mLf

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 :

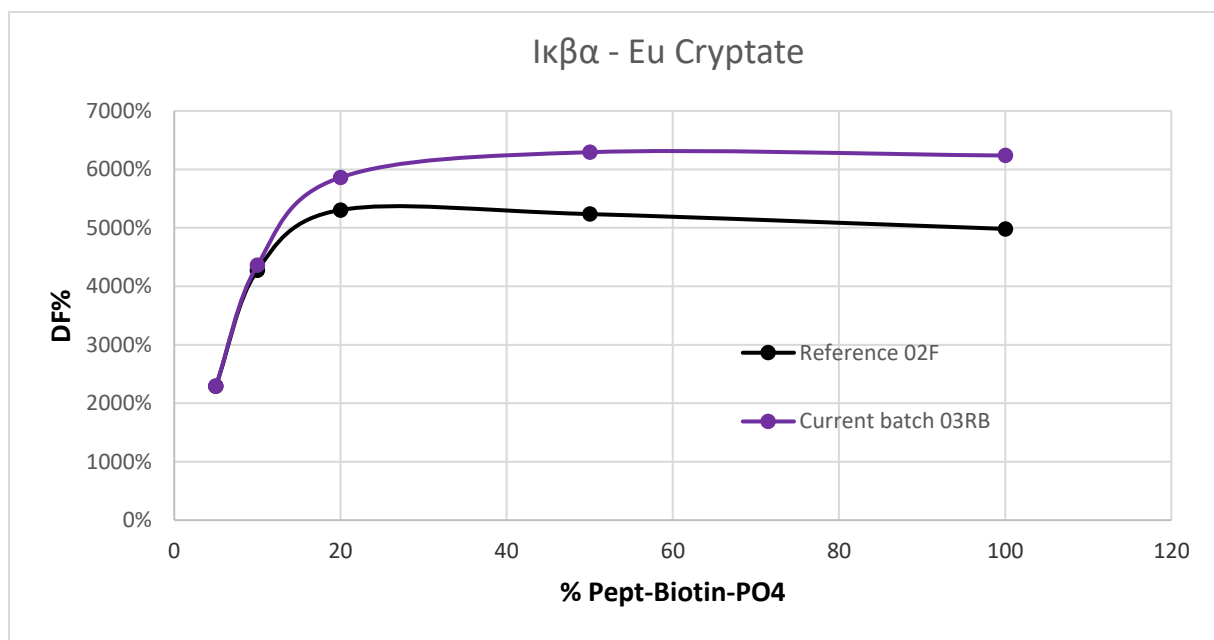
- 10 µL Peptide Biotin-PO4 and Peptide Biotin-non PO4
- 5 µL XL conjugate
- 5 µL K conjugate

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

## RESULTS

Peptide Biot-PO4 (n°839) %		Anti phospho IκBa -K Reference	Anti phospho IκBa-K CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
5	% Delta F % CV	2289 % 3.8 %	2294 % 2.7 %	0.2 %	11.6 %
10	% Delta F % CV	4272 % 0.2 %	4362 % 1.2 %	2.1 %	
20	% Delta F % CV	5305 % 0 %	5862 % 0.3 %	10.5 %	± 20 %
50	% Delta F % CV	5235 % 0.5 %	6294 % 0 %	20.2 %	
100	% Delta F % CV	4982 % 0.4 %	6237 % 0.1 %	25.2 %	

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



Anti phospho IκBa -K batch # 03RB

Accepted



Rejected



Operator's initials: LC

Date : 05/08/2005

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