

HTRF® Research Reagents

## FLAGM2-K Quality Control

Product number : 61FG2KLA Lot Number : 28RA

Manufacturing date : 06/03/2017 Document version : 1

### EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing Prot-N-Flag-Biot concentrations.

FLAGM2-K is lyophilized in PO4 buffer 100 mM pH7 ; 0.5% BSA ; Stabilizers.

All the reagents involved were diluted in 50 mM Hepes pH 8.0, 0.1% BSA, 0.4M KF.

### ASSAY FORMAT

FLAGM2-K + Prot-N-Flag-Biot + SA-XL

### REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
SA-XL	135	0.625 - 1.25 - 2.5 nMf
FLAGM2-K reference	25B	70 ng/mLf
Prot-N-Flag-Biot	CQ09	1.25 - 2.5 - 5 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
FLAGM2-K current batch	28RA	75 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

Reconstitute FLAGM2-K vial with 250 µL H2O (if packaged under 5,000 tests or 1,000 tests) or 1 mL H2O (if packaged under 20,000 tests).

Reagents were dispensed as follows:

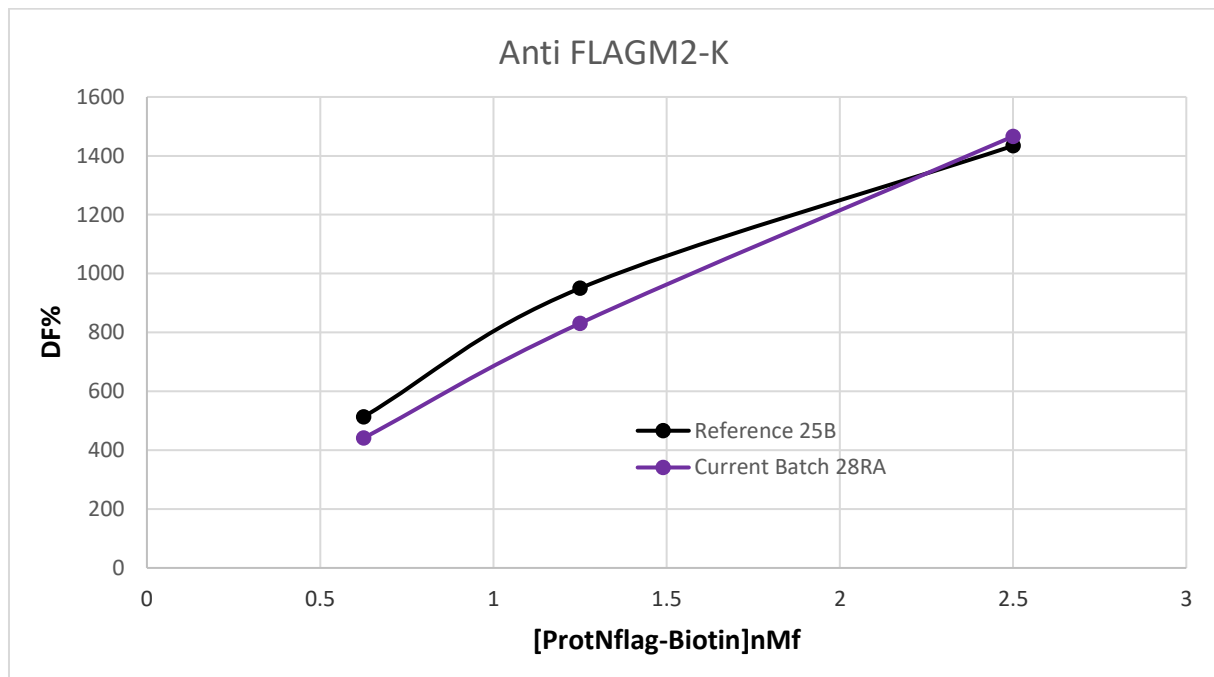
- 5 µL Prot-N-Flag-Biot (5 µL diluent for negative control)
- 10 µL XL conjugate
- 5 µL K conjugate

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

## RESULTS

[Prot-N-Flag-Biot] nMf		FLAGM2-K Reference	FLAGM2-K CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
1.25	% Delta F % CV	513% 2.0%	441% 1.0%	-14.0%	-8.1%
2.5	% Delta F % CV	950% 0%	831% 0.9%	-12.5%	± 20 %
5	% Delta F % CV	1434% 0%	1466% 1.0%	2.2%	

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



FLAGM2-K batch # 28RA

Accepted



Rejected



Operator's initials: MB

Date : 06/03/2017

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