

HTRF® Research Reagents

# Mab Anti Flag M2-K

# Quality Control

Product number : 61FG2KLA

Lot Number : 25C

Manufacturing date : 16/10/2024

Document version : 1

## EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing Prot-6HIS-Flag concentrations.

Mab Anti FlagM2-K is lyophilized in PO4 buffer 100 mM pH7; 0.5% BSA + Stabilizers.

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

## ASSAY FORMAT

Prot-6HIS-flag + 6HIS-XL + Mab Anti FlagM2-K

## REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
6HIS-XL	73A	1000 ng/mLf
FlagM2-K	29AR	75 ng/mlf
prot -6HIS- Flag	CQ11	0.156-0.62 - 2.5 -10 nMf
	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
FlagM2-K current batch	25C	70 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 Mab Anti FlagM2-Tb 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 :

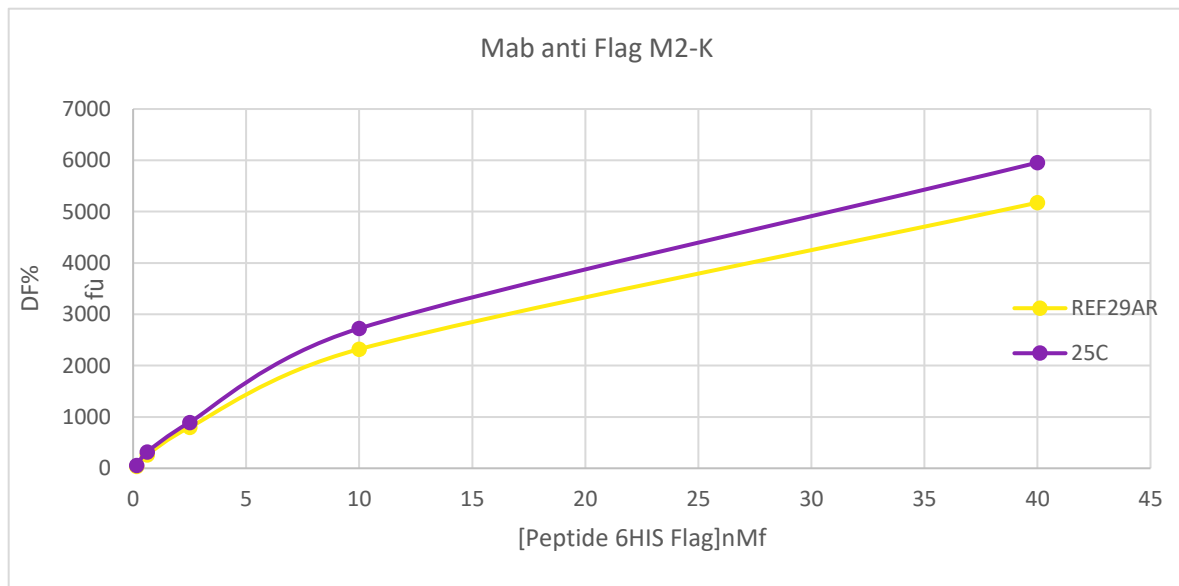
- 10 µL Prot-6HIS-Flag (10µL diluent for negative control)
- 5 µL XL conjugate
- 5 µL K conjugate

Incubation took place for Over Night at 4°C. The plate was read on Envision® under standard conditions.

## RESULTS

[Prot-6HIS -Flag] nMf		FlagM2-K Ref	FlagM2-K CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
0.156	% Delta F % CV	35 % 5.9%	55% 1.7 %		<b>16.8 %</b>
0.625	% Delta F % CV	262 % 0.8%	321 % 0.7%	22.5 %	
2.5	% Delta F % CV	798% 4.5%	895 % 2.1 %	12.2 %	± 20 %
10	% Delta F % CV	2321% 0.2%	2724% 1.1%	17.4 %	
40	% Delta F % CV	5175% 1.2%	5956% 0.8%	15.1%	

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



Flag M2-K batch # 25C

Accepted



Rejected



Operator's initials: ML

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