

HTRF® Research Reagents

## Anti FLAG-XL 665 Quality Control

Product number : 61FGBXLB Lot Number : 01A

Manufacturing date : 18/09/2024 Document version : 1

### EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing antigen concentrations (Biot-6HIS-FLAG).

anti Flag-XL is lyophilized in PO4 buffer 100 mM pH7 ; 0.5% BSA ; Stabilizers.

All the reagents involved were diluted in HTRF PPI-Terbium detection buffer (#61DB10RDF).

### ASSAY FORMAT

anti-6HIS-Tb GOLD + Biot-6HIS-FLAG + anti-FLAG-XL665

### REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
Anti-6HIS-Tb GOLD	19A	0.3 nMf
Anti-FLAG-XL665 référence	-	6.67 nMf
Biot-6HIS-FLAG	01	10, 5, 2.5, 1.25, 0.62, 0.31 et 0.15nM nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti-FLAG-XL665 current batch	01A	6.67 nMf

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

### EXPERIMENTAL PROCEDURE

Reconstitute anti Flag-XL665 vial (20,000 tests) with 1 mL H2O.

After reconstitution:

-Concentration (total protein) : 0.4 mg/mL

Reagents were dispensed as follows :

- 10 µL Biot-6HIS-FLAG (10 µL diluent for negative control)
- 5 µL XL conjugate
- 5 µL K conjugate Tb

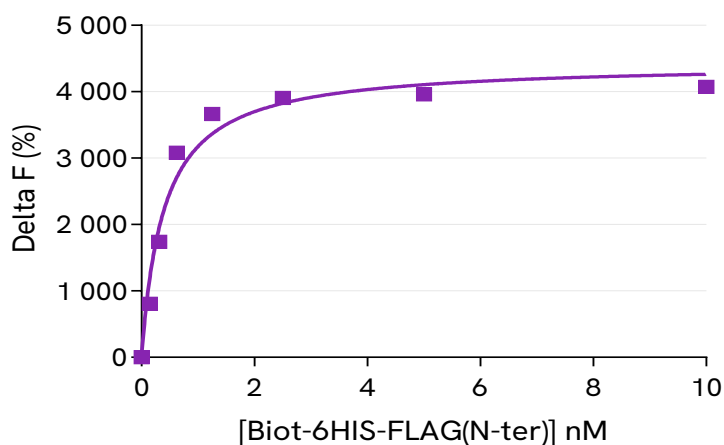
Incubation took place for 120 min at room temperature. The plate was read on Envision under standard conditions.

## RESULTS

[Biot-6HIS-FLAG] nMf		anti Flag-XL Reference	anti Flag-XL CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
0.15	% Delta F % CV	-	807 % 3.7 %	-	
0.31	% Delta F % CV	-	1739 % 3.2 %	-	± 20 %
0.62	% Delta F % CV	-	3080 % 1 %	-	
1.25	% Delta %cv	-	3665% 0.1%	-	
2.5	% Delta %cv	-	3905 % 0.6%	-	
5	% Delta %cv	-	3962% 1.2%	-	
10	% Delta %cv	-	4075 % 1.1%	-	

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

### Toolbox anti FLAG-XL665



	Kd
M2	0.3965

**Kd accepted range**  
**0,2416 to 0,7046 nM**

anti Flag-XL 665 batch # 01A

Accepted



Rejected



Operator's initials: FL

Date: 2024-12-05

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