

HTRF® Research Reagents

## Anti-6HIS-XL665 Quality Control

Product number : 61HISXLF Lot Number : 70RD

Manufacturing date : 29/09/2021 Document version : 1

### EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing Anti-6HIS-XL665 concentrations.

Anti-6HIS-XL665 is lyophilized in PO4 buffer 100 mM pH7 ; 0.1% BSA ; Stabilizers.

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

### ASSAY FORMAT

Pept-6HIS-Biotin + Streptavidin-Cryptate + Anti-6HIS-XL665

### REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
Streptavidin-Cryptate	34A	1.8 nMf
Anti-6HIS-XL665 reference	66B	250 à 1 000 ng/mLf
Pept-6HIS-Biotin	CQ02	5 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti-6HIS-XL665 current batch	70RD	250 à 1 000 ng/mLf

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

### EXPERIMENTAL PROCEDURE

Reconstitute Anti-6HIS-XL665 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 :

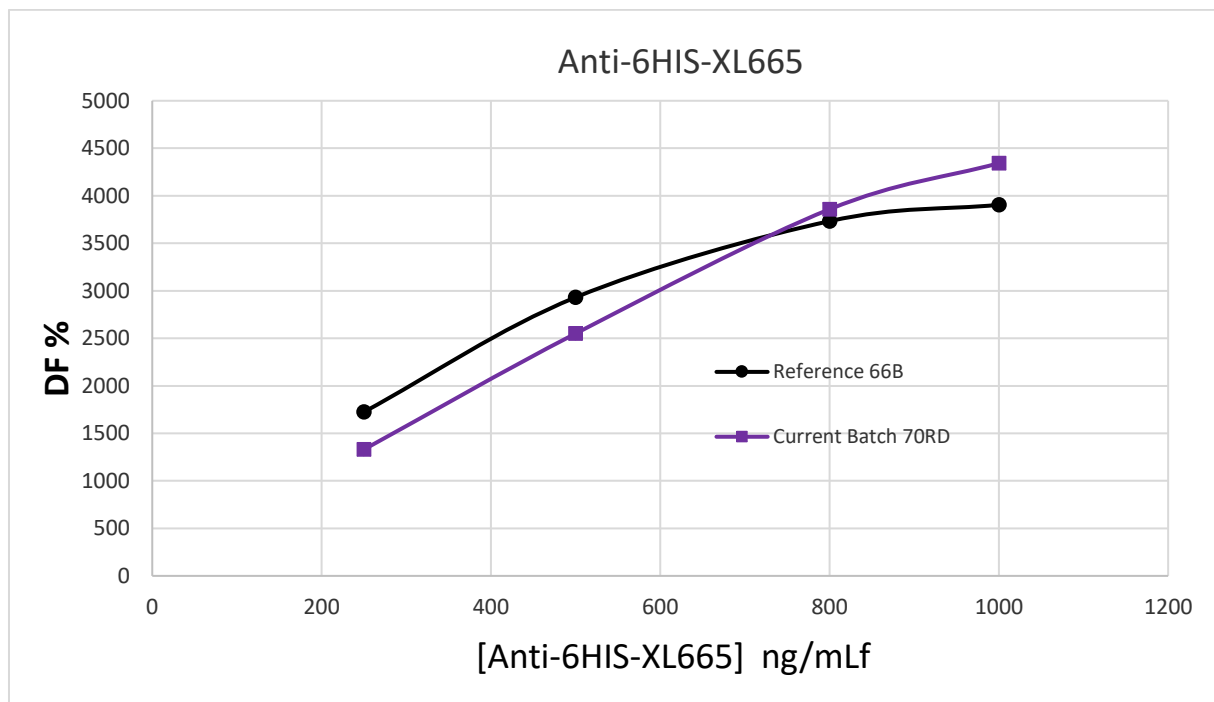
- 5 µL Pept-6HIS-Biotin (5 µL diluent for negative control)
- 10 µL XL conjugate
- 5 µL K conjugate

Incubation took place for 150 min at 2-8°C. The plate was read on Phera Star ® under standard conditions.

## RESULTS

[Anti-6HIS-XL665] ng/mLf		Anti-6HIS-XL665 Reference	Anti-6HIS-XL665 CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
250	% Delta F % CV	1725 % 1.7 %	1330 % 0.7 %	-22.9 %	-5.3 %
500	% Delta F % CV	2931 % 1.7 %	2550 % 1.0 %	-13.0 %	
800	% Delta F % CV	3733 % 1.8 %	3858 % 1.3 %	3.3 %	± 20 %
1000	% Delta F % CV	3905 % 1.8 %	4343 % 3.9 %	11.2 %	

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



Anti-6HIS-XL665 batch # 70RD

Accepted

Rejected

Operator's initials: AL

Date: 29/09/2021

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