

HTRF® Research Reagents

Mab Anti-6HIS-Eu cryptate Gold Quality Control

Product number : 61HI2KLB Lot Number : 18A

Manufacturing date : 18/01/2024 Document version : 1

EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing peptide concentrations (Pept-6HIS-Biotin).

MAB Anti-6HIS-Eu cryptate Gold 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

MAB Anti-6HIS-Eu cryptate Gold + Pept-6HIS-biot. + Streptavidin-XL665

REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
Streptavidin-XL665	181A	2.4 nMf à 0.15 nMf
MAB Anti-6HIS-Eu cryptate Gold reference	17A	0.5 nMf
Pept-6HIS-Biot.	CQ03	9.6 nMf à 0.6 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
MAB Anti-6HIS-Eu cryptate Gold current Batch	18A	0.5 nMf

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

If applicable, streptavidin / biotin ratio was set up at : 1/4

EXPERIMENTAL PROCEDURE

Reconstitute MAB Anti-6HIS-Eu cryptate Gold 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 :

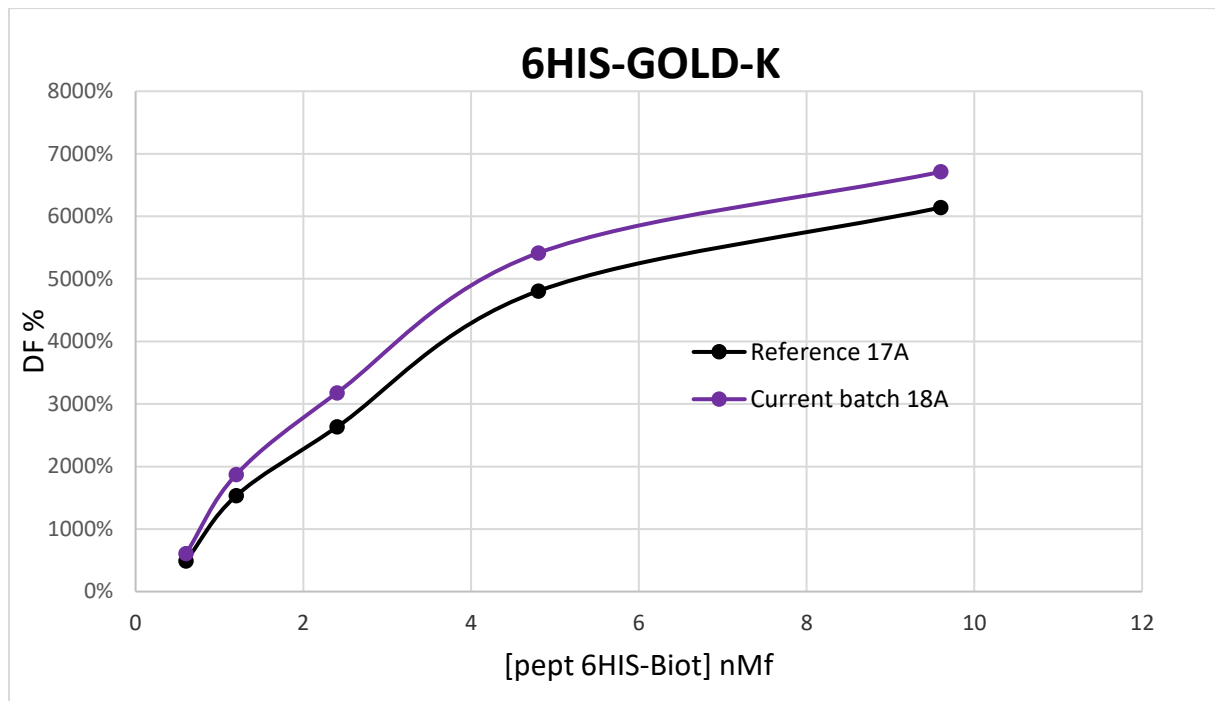
- 10 µL Pept 6HIS-Biot
- 5 µL XL conjugate
- 5 µL Eu conjugate

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

RESULTS

[Pept-6HIS-Biot] nM _f		Anti-6HIS-Eu cryptate Gold Reference	Anti-6HIS-Eu cryptate Gold Current batch	Variation / Ref	Mean Acceptable variation
0.6	% Delta F	488%	603%	23.7%	
	% CV	0.6%	2.0%		
1.2	% Delta F	1532%	1870%	22.1%	
	% CV	0.5%	1.0%		
2.4	% Delta F	2630%	3176%	20.8%	
	% CV	1.8%	1.0%		
4.8	% Delta F	4805%	5413%	12.6%	
	% CV	2.6%	1.4%		
9.6	% Delta F	6140%	6710%	9.3%	± 20 %
	% CV	1.1%	1.3%		

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



Anti-6HIS-Eu cryptate Gold # 18A

Accepted

Rejected

Operator's initials: FMV

Date : 31/01/2024

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.

