

HTRF® Research Reagents

## Anti-6His-D2 Quality Control

Product number : 61HISDLA Lot Number : 39

Manufacturing date : 08/05/2023 Document version : 1

### EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently used, at increasing mab concentrations (Anti-6HIS-d2).

Anti-6HIS-d2 is lyophilized in PO4 buffer 100 mM pH 7 ; 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-d2

### REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
Streptavidin-Cryptate	37A	110 ng/ml <sub>f</sub>
Anti-6HIS-d2	34A	125, 250, 500, 800 ng/ml <sub>f</sub>
Pept-6HIS-Biotin	CQ03	5 nM <sub>f</sub>

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti-6HIS-d2	39	125, 250, 500, 800 ng/ml <sub>f</sub>

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

### EXPERIMENTAL PROCEDURE

Reconstitute Anti-6HIS-d2 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 d2 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-d2] nMf		anti -HIS-d2 Ref	anti 6HIS-d2 CURRENT BATCH	VARIATION	ACCEPTABLE VARIATION
125	% Delta F % CV	850 % 0.8%	899 % 6%	5.8%	5.6 %
250	% Delta F % CV	1351 % 1.7%	1481 % 10 %	9.7 %	
500	% Delta F % CV	1688 % 4.5 %	1762 % 4.4 %	4.4 %	Acceptable variation
800	% Delta F % CV	1766 % 6.7 %	1813 % 2.7 %	2.7 %	+/- 20 %

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

Anti 6HIS-d2 batch # 39A

Accepted

Rejected

Operator's initials:

FMV

20/08/2024

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.

revvity