

HTRF® Research Reagents

## Anti HA-d2 Quality Control

Product number : 610HADAA

Lot Number : 14BR

Manufacturing date : 19/09/2023

Document version : 1

### EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing anti HA-d2 concentrations.

Anti HA-d2 is frozen in PO4 buffer 100 mM pH7; 0.1% BSA .

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

### ASSAY FORMAT

Pept-HA-DNP + Anti HA-d2 + Anti DNP -Cryptate

### REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
Anti DNP Cryptate	15A	200 ng/mLf
Anti HA-d2 Ref	14RA	0.0156 - 1 nMf
Peptide HA-DNP	CQ02	1 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti HA-d2 S current batch	14BR	0.0156 - 1 nMf

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

### EXPERIMENTAL PROCEDURE

Reagents were dispensed as follows :

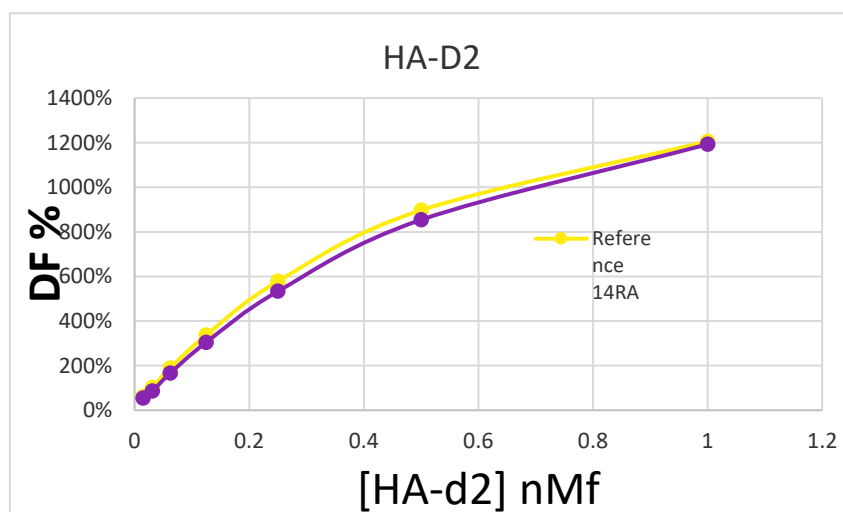
- 5 µL Peptide HA-DNP (5 µL diluent for negative control)
- 10 µL d2 conjugate (anti HA-d2)
- 5 µL K conjugate ( Anti DNP-Cryptate)

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

## RESULTS

[Peptide HA-DNP] nMf		HA-d2 Ref	HA-d2 CURRENT BATCH	VARIATION	ACCEPTABLE VARIATION
0.0156	% Delta F % CV	62 % 7.7 %	54 % 7.2%	-11.5 %	-9.01 %
0.0313	% Delta F % CV	102 % 1.7 %	86 % 0.9%	-15.8 %	
0.625	% Delta F % CV	189 % 1.3 %	166 % 1.6 %	-12.5 %	± 20 %
0.125	% Delta F % CV	337% 0.9 %	304% 0.5 %	-9.6 %	
0.25	% Delta F % CV	578 % 0.6 %	533 % 3.0 %	-7.7%	
0.5	% Delta F % CV	894 % 7.0 %	854 % 1.2 %	-4.8%	
1	% Delta F % CV	1206 % 0.2%	1192 % 1.1 %	-1.1 %	

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



HA-d2 batch # 14BR

Accepted



Rejected



Operator's initials: ML

Date : 22/09/2023

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