

HTRF® Research Reagents

## ANTI hFc-d2 Quality Control

Product number : 61HFCDAF Lot Number : 11A

Manufacturing date : 08/11/2024 Document version : 1

### EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing anti body concentrations (Ac Human-Biotin).

Anti hFc-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

Ac Human Biotin + Anti hFc -d2 + SA-K

### REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
SA-K	37 A	110 ng/mLf
Anti hFc-d2 Ref	10A	1000 nMf
Ac Human-Biotin	CQ01	0.187 - 1.5 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti hFc-d2 current batch	11A	1000 nMf

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

### EXPERIMENTAL PROCEDURE

Reagents were dispensed as follows :

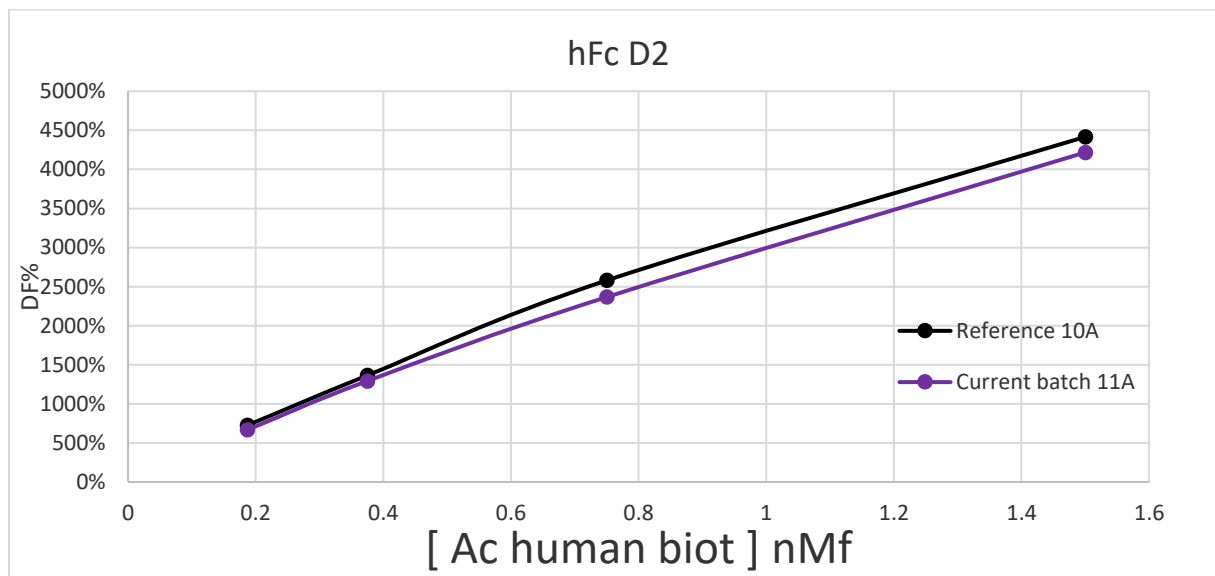
- 5 µL Ac Human-Biotin
- 10 µL d2 conjugate
- 5 µL K conjugate

Incubation took place for 180 min at room temperature. The plate was read on Phera Star ® under standard conditions.

## RESULTS

[Ac Human-Biotin] nMf		Anti hFc-d2 Ref	Anti hFc-d2 CURRENT BATCH	VARIATION	ACCEPTABLE VARIATION
0.187	% Delta F % CV	725 % 3.6 %	667% 4.4 %	-8.1 %	-6.5%
0.375	% Delta F % CV	1363 % 1.0 %	1292% 1.3 %	-5.2%	
0.75	% Delta F % CV	2579 % 3.8 %	2366 % 3.1 %	-8.3 %	± 20 %
1.5	% Delta F % CV	4416 % 2.9 %	4217 % 2.7 %	-4.5 %	

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



Anti hFc-d2 batch # 11A

Accepted

Rejected

Operator's initials: ML

Date : 12/11/2024

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