

HTRF® Research Reagents

Anti hFc-XL Quality Control

Product number : 61HFCXLA Lot Number : 14RA

Manufacturing date : 15/09/2022 Document version : 1

EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing Ab Human-Biotin concentrations.

Anti hFc-XL 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

Ab Human-Biotin + Anti hFc-XL + SA-K

REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
SA-K	34A	110 ng/mLf
Anti hFc-XL reference	13A	1000 ng/mLf
Ab Human-Biotin	CQ01	0.187 nMf – 1.5 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti hFc-XL current batch	14RA	1000 ng/mLf

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

EXPERIMENTAL PROCEDURE

Reconstitute Anti hFc-XL 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 :

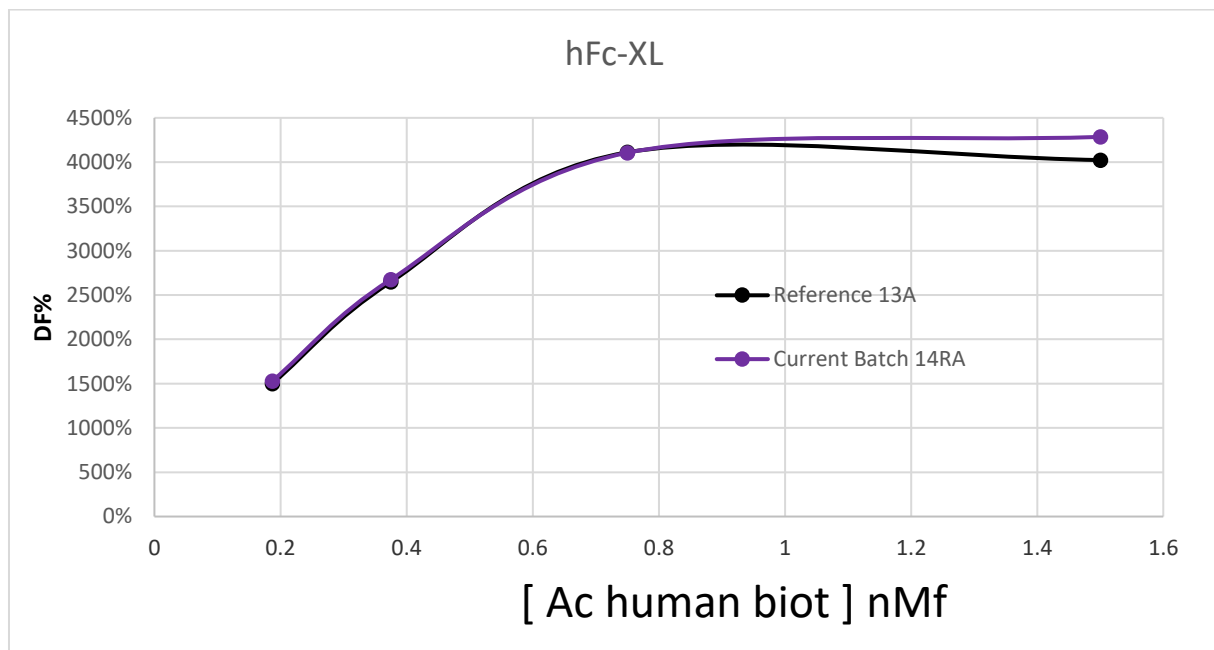
- 5 µL Ab Human-Biotin (5 µL diluent for negative control)
- 10 µL XL conjugate
- 5 µL K conjugate

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

RESULTS

[Ab Human-Biotin] nMf		Anti hFc-XL Reference	Anti hFc-XL CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
0.187	% Delta F % CV	1501 % 2.9 %	1530 % 1.1 %	2.0 %	2.3 %
0.375	% Delta F % CV	2647 % 3.4 %	2675 % 0.7 %	1.0 %	
0.75	% Delta F % CV	4110 % 1.5 %	4106 % 0.9 %	-0.1 %	± 20 %
1.5	% Delta F % CV	4023 % 1.6 %	4285 % 1.3 %	6.5 %	

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



Anti hFc-XL batch # 14RA

Accepted



Rejected



Operator's initials: MB

Date : 16/09/2022

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