

HTRF® Research Reagents

Anti-Rabbit-XL Quality Control

Product number : 61PARXLB Lot Number : 13RA

Manufacturing date : 2018-03-19 Document version : 1

EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing Anti-Rabbit-XL concentrations.

Anti-Rabbit-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

SA-K + Ab Rabbit-Biotin + Anti-Rabbit-XL

REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
SA-K	CQ02	2.5 nMf
Anti Rabbit-XL reference	12A	1.67 - 3.34 - 6.67 nMf
Ab Rabbit-Biotin	CQ01	1.67 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Anti-Rabbit-XL current batch	13RA	1.67 - 3.34 - 6.67 nMf

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

EXPERIMENTAL PROCEDURE

Reconstitute Anti-Rabbit-XL vial with 250 µL H2O (if packaged under 5 000 tests) or 1 mL H2O (if packaged under 20 000 tests).

Reagents were dispensed as follows :

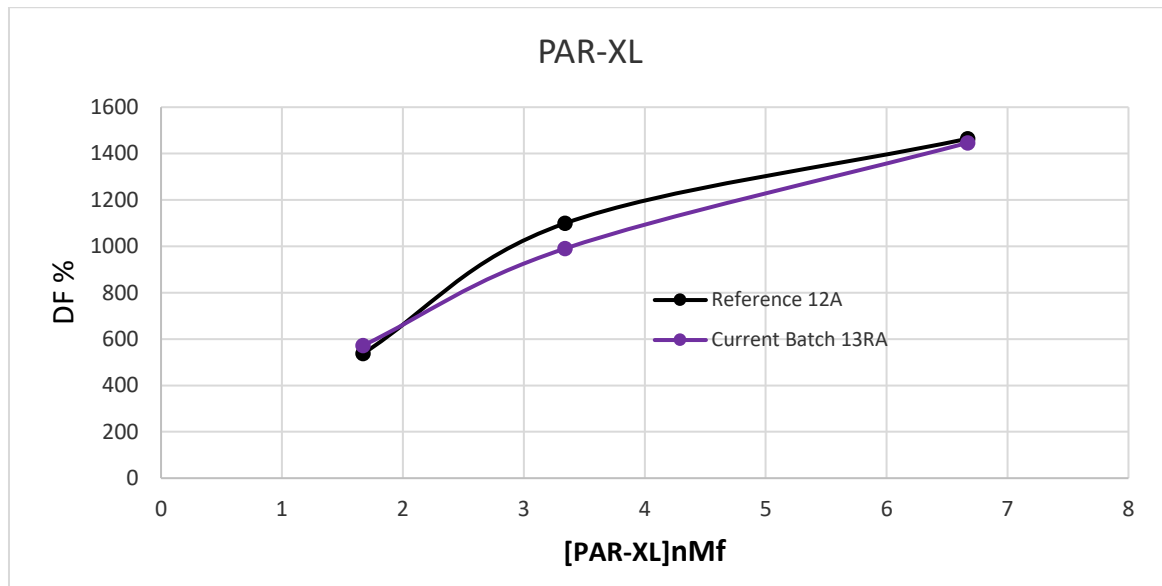
- 5 µL PPI - Europium detection buffer
- 5 µL Ab Rabbit-Biotin
- 5 µL XL conjugate
- 5 µL K conjugate

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

RESULTS

[Anti-Rabbit-XL] nM _f		Anti-Rabbit-XL Reference	Anti-Rabbit-XL CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
1.67	% Delta F % CV	537 % 1.1 %	572 % 1.6 %	6.5 %	-1.5 %
3.34	% Delta F % CV	1 099 % 0.3 %	990 % 1.2 %	-9.9 %	
6.67	% Delta F % CV	1 464 % 8.6 %	1 446 % 7.7 %	-1.2 %	± 20 %

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



Anti-Rabbit-XL batch # 13RA

Accepted

Rejected

Operator's initials: FTV

Date: 16/03/2018

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