

HTRF® Research Reagents

## Protein A-XL665 Quality Control

Product number : 61PRAXLB Lot Number : 19A

Manufacturing date : 04/11/2024 Document version : 1

### EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently used, at increasing protein A-XL concentrations.

Protein A-XL is lyophilized in PO<sub>4</sub> buffer 100 mM pH7, 0.1% BSA ; Stabilizers.

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

### ASSAY FORMAT

Protein A-XL + anti GST-Cryptate (or Streptavidin-Cryptate)

### REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
Streptavidin-Cryptate	29A	1.5 nM <sub>f</sub>
Protein A-XL	17A	0.125 nM <sub>f</sub> - 2 nM <sub>f</sub>
Anti GST-Cryptate	107A	0.83 nM <sub>f</sub>

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
Protein A-XL	19A	0.125 nM <sub>f</sub> - 2 nM <sub>f</sub>

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

### EXPERIMENTAL PROCEDURE

Protein A-XL vial was reconstituted with 250 µl H<sub>2</sub>O (if packaged under 5,000 tests) or 1 ml H<sub>2</sub>O (if packaged under 20,000 tests).

Reagents were dispensed as follows :

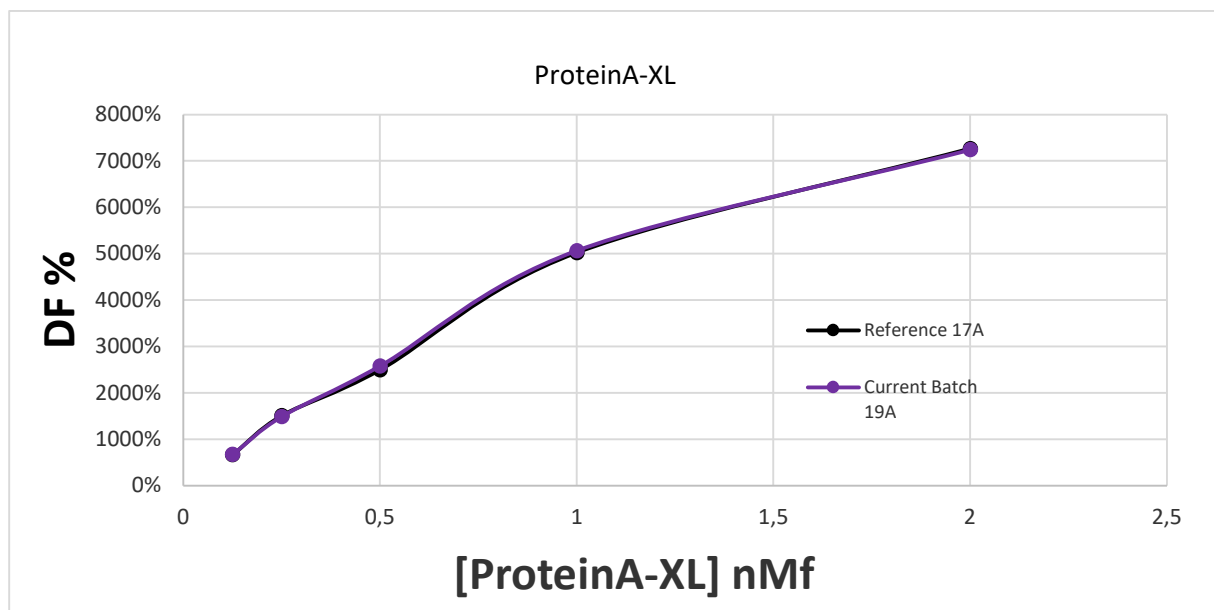
- 10 µl Cryptate conjugate
- 10 µl XL conjugate

Incubation took place for 60 min at room temperature. The plate was read on Phera Star Plus® under standard conditions.

## RESULTS

[Protéine A-XL] nM <sub>f</sub>		Protéine A-XL Réf	Protéine A-XL 20000 pts Current batch	VARIATION	MEAN ACCEPTABLE VARIATION
0.125	% Delta F % CV	666% 0.7 %	675% 0.2 %	1.3%	<b>0.72%</b>
0.25	% Delta F % CV	1512 % 1.5 %	1492% 3.1 %	-1.3%	
0.5	% Delta F % CV	2500 % 2.5 %	2580 % 4.0%	3.2%	± 20 %
1	% Delta F % CV	5026 % 0.8 %	5065% 1.2 %	0.8%	
2	% Delta F % CV	7268 % 1.4 %	7241 % 0.8%	-0.4%	± 20 %

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



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

Protein A-XL batch # 19A

Accepted

Rejected

Operator's initials: MP

Date : 12/11/2024