

HTRF® Research Reagents

PolyGT-BIOTIN Quality Control

Product number : 61GT0BLA Lot Number : 32RA

Manufacturing date : 07/04/2020 Document version : 1

EXPERIMENTAL CONDITION

The batch to be controlled was compared to the reference batch currently in use, at increasing PolyGT-Biotin concentrations and SA-XL concentrations.

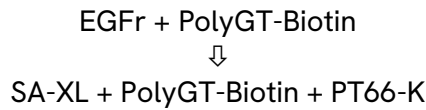
Phosphorylation of the peptide was carried using EGFr in Enzymatic Buffer (62EZBFDC) 1X + enzymatic cofactor for 20 min.

Revelation by SA-XL and PT66-K was carried out for 20 min.

PolyGT-Biotin is lyophilized in 20 mM Hepes pH 7.4 ; 1% BSA ; Stabilizers.

All revelation reagents involved were diluted in 50 mM Hepes pH 7, 0.1% BSA, 0.5M KF.

ASSAY FORMAT



REAGENTS

	REFERENCE LOT NUMBER	ASSAY CONCENTRATION
SA-XL	158A	2.5 nMf - 100 nMf
PT66-K	52B	125 ng/mLf
PolyGT-Biotin reference	31A	2.5 nMf - 100 nMf

	CURRENT BATCH LOT NUMBER	ASSAY CONCENTRATION
PolyGT-Biotin current batch	32RA	2.5 nMf - 100 nMf

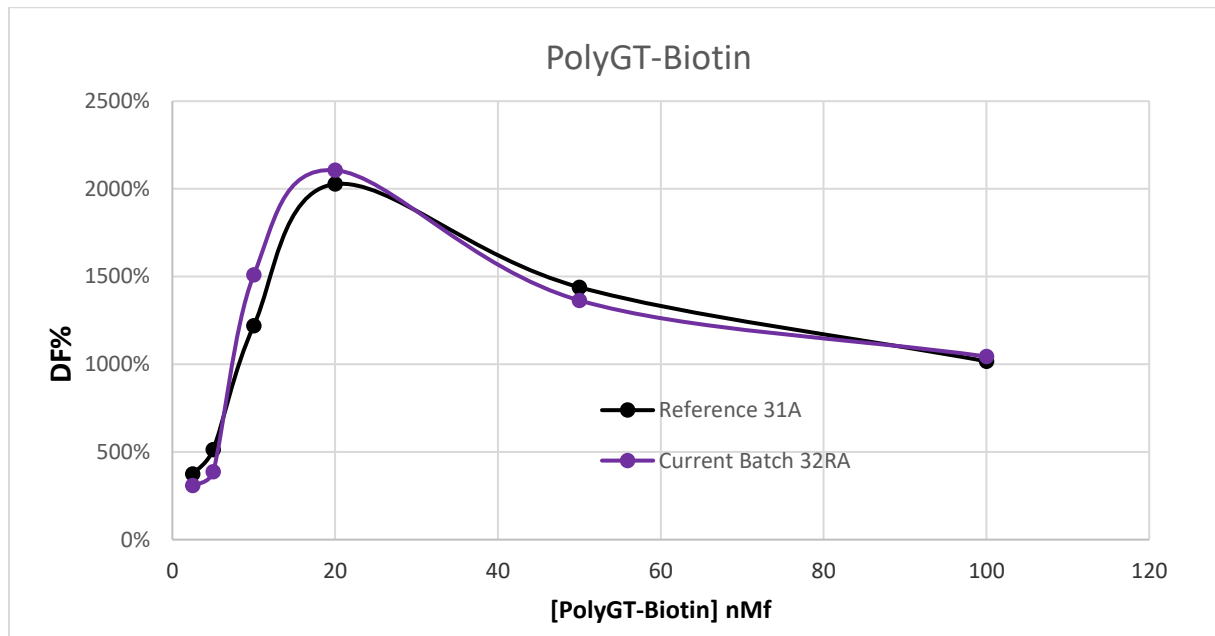
Other reagents were the same as those used for the reference test.

The plate was read on Phera Star ® under standard conditions.

RESULTS

[PolyGT-Biotin] nMf		PolyGT-Biotin Reference	PolyGT-Biotin CURRENT BATCH	VARIATION	MEAN ACCEPTABLE VARIATION
2.5	% Delta F % CV	374 % 19.1 %	309 % 1.4 %	-2.2 %	
5	% Delta F % CV	513 % 6.1 %	386 % 0 %	-24.9 %	-0.3 %
10	% Delta F % CV	1220 % 0 %	1509 % 1.4 %	23.7 %	
20	% Delta F % CV	2027 % 14.4 %	2106 % 4.9 %	3.9 %	
50	% Delta F % CV	1439 % 22.5 %	1364 % 1.8 %	-5.2 %	± 20 %
100	% Delta F % CV	1017 % 0.9 %	1044 % 4.3 %	2.7 %	

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



PolyGT-Biotin batch # 32RA

Accepted

Rejected

Operator's initials: MB

Date : 07/04/2020

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.

