

HTRF® Research Reagents

Anti C-MYC-Tb Quality Control

Product number : 61MYCTAF Lot Number : 04RB

Manufacturing date : 08/02/2018 Document version : 1

EXPERIMENTAL CONDITION

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

Anti C-MYC-Tb is frozen in PO4 buffer 100 mM pH7 ; 0.1% BSA .

All the reagents involved were diluted in PPI Terbium detection buffer (#61DB10RDF).

ASSAY FORMAT

Anti C-MYC-Tb + Peptide C-MYC-Biotin + SA-XL

REAGENTS

| | REFERENCE LOT NUMBER | ASSAY CONCENTRATION |
|-------------------------|----------------------|--------------------------|
| SA-XL | 135A | 0.312-0.625-1.25-2.5 nMf |
| Anti C-MYC-Tb reference | 03B | 0.5 nMf |
| Peptide C-MYC-Biotin | CQ01 | 0.312-0.625-1.25-2.5 nMf |

| | CURRENT BATCH LOT NUMBER | ASSAY CONCENTRATION |
|-----------------------------|--------------------------|---------------------|
| Anti C-MYC-Tb current batch | 04RB | 0.5 nMf |

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

If applicable, Biotin/Streptavidin ratio was set up at: 1/1

EXPERIMENTAL PROCEDURE

Reagents were dispensed as follows :

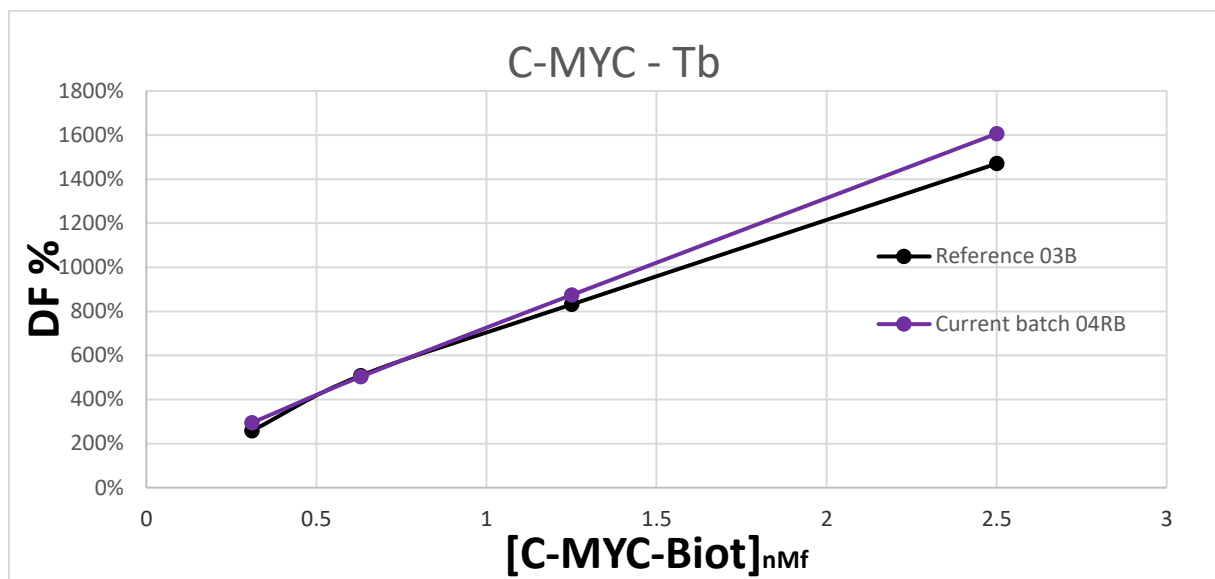
- 5 µL Peptide C-MYC-Biotin (5 µL diluent for negative control)
- 10 µL XL conjugate
- 5 µL Tb conjugate

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

RESULTS

| [SA-XL] nMf | | Anti C-MYC-Tb Reference | Anti C-MYC-Tb Current batch | VARIATION | ACCEPTABLE VARIATION |
|----------------|-------------------|----------------------------|--------------------------------|-----------|-------------------------|
| 0.3125 | % Delta F % cv | 258 % 1.9 % | 294 % 0.6 % | 14.0 % | 6.9 % |
| 0.625 | % Delta F % cv | 508 % 0 % | 503 % 2.4 % | -1.0 % | |
| 1.25 | % Delta F % cv | 831 % 2.6 % | 874 % 1.2 % | 5.2 % | +/- 20% |
| 2.5 | % Delta F % cv | 1470 % 0.4 % | 1606 % 2.9 % | 9.2 % | |

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



Anti-C-MYC-Tb batch # 04RB

Accepted



Rejected



Operator's initials: MB

Date: 08/02/2018

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