

## HTRF® Research Reagents

**HTRF P53 total kit - 96 pts**

Product Number : 64T53PET Lot Number : 04F

Expiration date : 2024-09-09 Document version : 1

| Reagents :                                       |  | Lot n° |
|--|--|--------|
| PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND           |  | 2203   |
| HTRF P53 total kit - Ctrl lysate                 |  | 04RA   |
| HTRF P53 total kit - 96 pts d2 antibody          |  | 04RB   |
| HTRF P53 total kit - 96 pts Eu Cryptate antibody |  | 04RB   |
| HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL           |  | 35AR   |
| HTRF P-T prot. - Block. reag.(100X) 0.3 mL       |  | 28A    |
| HTRF P-T prot. - Detect. Buf. 0.5 mL             |  | 23RB   |

**Results :**

 Incubation : Overnight at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Pherastar Plus  
 Fitting model : /

**Assay characteristics :**

|                | Found value | Specifications Rev 01 |
|----------------|-------------|-----------------------|
| Negative       | 428         | /                     |
| Control Lysate | 3374        | /                     |
| S/B            | 7,9         | >5                    |
|                | /           | /                     |

**HTRF P53 total kit - 96 pts - Lot n° : 04F**
 Accepted

 Rejected

Quality control date : 2022-03-01 Supervisor's initials : AC

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.