

HTRF® Research Reagents

## HTRF P38 p-T180/Y182 kit - 500 pts

Product Number : 64P38PEG Lot Number : 07C

Expiration date : 07/12/2025 Document version : 1

| Reagents :  | Lot n° |
|---|--------|
| HTRF P38 p-T180/Y182 kit - 500 pts d2 antibody          | 07RA   |
| HTRF P38 p-T180/Y182 kit - 500 pts Eu Cryptate antibody | 07RA   |
| HTRF P38 p-T180/Y182 kit - Ctrl lysate                  | 12RA   |
| HTRF P38 p-T180/Y182 kit - modop                        | 7      |
| HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL                  | 33A    |
| HTRF P-T prot. - Block. reag.(100X) 0.3 mL              | 28A    |
| HTRF P-T prot. - Detect. Buf. 2 mL                      | 32A    |

**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       | 563         | /                     |
| Control Lysate | 6554        | /                     |
| S/B            | 11,6        | >5                    |
|                | /           | /                     |

HTRF P38 p-T180/Y182 kit - 500 pts - Lot n° : 07C

 Accepted

 Rejected

|                                   |                            |
|-----------------------------------|----------------------------|
| Quality control date : 2023-06-01 | 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.