

HTRF® Research Reagents

HTRF KRAS WT/SOS1 PPI kit - 500 pts

Product Number : 64KRASWTPEG Lot Number : 02D

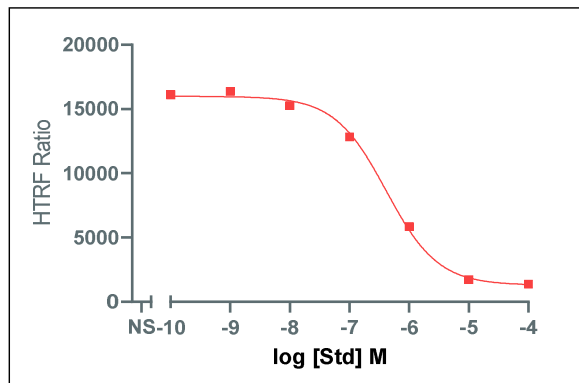
Expiration date : 2025-02-24 Document version : 1

| Reagents : | Lot n° |
|--|--------|
| HTRF KRAS WT/SOS1 PPI kit - 500 pts Tag2 Human SOS1 | 02RA |
| HTRF KRAS/SOS1 PPI kit - Std | 03RA |
| HTRF KRAS WT/SOS1 PPI kit - 500 pts GTP | 03RA |
| HTRF KRAS WT/SOS1 PPI kit - 500 pts Anti Tag2 Tb Cryptate antibody | 02RA |
| HTRF KRAS WT/SOS1 PPI kit - 500 pts Tag1 Human KRAS WT | 02RA |
| HTRF KRAS WT/SOS1 PPI kit - 500 pts Anti-tag1 XL665 antibody | 02RA |
| HTRF KRAS WT/SOS1 PPI kit - modop | 2 |
| HTRF Bind. Domain - Detect. Buf. 2- 20 mL | 11RA |

Results :

Incubation : 2h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : log(inhibitor) vs.response-Variable slope (four parameters)

| STD | Conc. (µMf) | RATIO |
|-------|-------------|--------|
| Std 0 | 0 | 16 106 |
| Std 1 | 0.001 | 16 353 |
| Std 2 | 0.01 | 15 280 |
| Std 3 | 0.1 | 12 844 |
| Std 4 | 1 | 5 861 |
| Std 5 | 10 | 1 722 |
| Std 6 | 100 | 1 371 |
| Std 7 | 1000 | 1 290 |



Assay characteristics :

| | Found value | Specifications Rev 01 |
|-------|-------------|-----------------------|
| S/B | 15 | 11-15 |
| IC 50 | 412 | 300-600 nM |

HTRF KRAS WT/SOS1 PPI kit - 500 pts - Lot n° : 02D

Accepted

Rejected

| | |
|-----------------------------------|----------------------------|
| Quality control date : 2023-10-17 | Supervisor's initials : FL |
|-----------------------------------|----------------------------|

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.