

HTRF® Research Reagents

HTRF IP-One Gq kit - 1K pts

Product Number : 62IPAPEB Lot Number : 24B

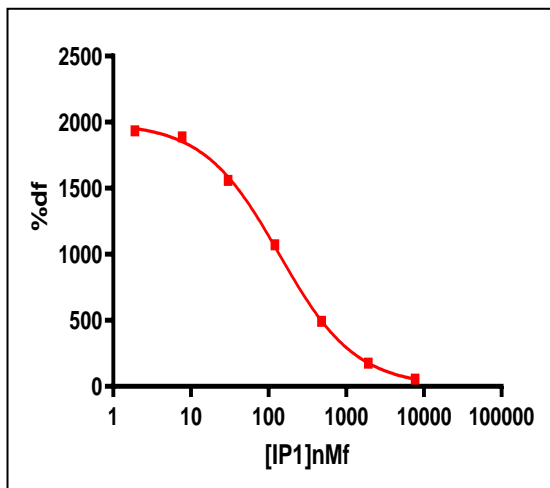
Expiration date : 2026-08-02 Document version : 1

| Reagents : | Lot n° |
|-------------------------------------|--------|
| HTRF Lysis & Detect. Buf. 5 - 13 mL | 08A |
| HTRF IP-One Gq kit - 1K pts Tb | 24A |
| HTRF IP-One Gq kit - 1K pts d2 | 24A |
| HTRF IP-One Gq kit - Std | 16AR |
| HTRF Stimulation Buf. 2 (5X) - 8 mL | 39A |
| / | / |
| / | / |

Results :

Incubation : 1h at RT
 Plate : 384 wells regular white plate 30 µl
 Reader : Pherastar Plus
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

| IP-One | Conc. (nMf) | DF % |
|--------|-------------|-------|
| Std 0 | 0 | 1885% |
| Std 1 | 1,9 | 1934% |
| Std 2 | 7,7 | 1889% |
| Std 3 | 30,1 | 1560% |
| Std 4 | 120,4 | 1071% |
| Std 5 | 481,6 | 492% |
| Std 6 | 1925 | 177% |
| Std 7 | 7700 | 56% |



Assay characteristics :

| | Found value | Specifications Rev 01 |
|----------|-------------|-----------------------|
| DF max % | 1885% | 1800 – 2600 % |
| EC 50 | 138,1 | 100 - 140 nMf |
| S/B | / | / |
| | / | / |

HTRF IP-One Gq kit - 1K pts - Lot n° : 24B

Accepted

Rejected

Quality control date : 2024-12-23 Supervisor's initials : MP

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.