

HTRF® Research Reagents

HTRF cAMP Gs HiRange kit - 100K pts

Product Number : 62AM6PEJ Lot Number : 21E

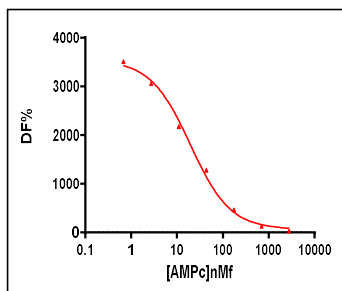
Expiration date : 2026-03-31 Document version : 1

| Reagents : | Lot n° |
|---------------------------------------|--------|
| HTRF Lysis & Detect. Buf. 2 - 200 mL | 31A |
| HTRF Stimulation Buf. 1 (5X) - 100 mL | 36A |
| HTRF cAMP Gs HiRange kit - Std | 19A |
| HTRF cAMP Gs HiRange kit - 20K pts Eu | 21A |
| HTRF cAMP Gs HiRange kit - 20K pts d2 | 21A |

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL (PROXIPLATE)
 Reader : Pherastar Plus
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

| cAMP | Conc. (nMf) | DF (%) |
|-------|-------------|--------|
| Std 0 | 0 | 3758 |
| Std 1 | 0,68 | 3509 |
| Std 2 | 2,73 | 3063 |
| Std 3 | 10,94 | 2167 |
| Std 4 | 43,75 | 1278 |
| Std 5 | 175 | 458 |
| Std 6 | 700 | 124 |
| Std 7 | 2800 | 29 |



Assay characteristics :

| | Found value | Specifications Rev 01 |
|---------|-------------|-----------------------|
| DF% max | 3758% | 2900 % - 4150% |
| EC 50 | 17,42 | 10 nMf - 22 nMf |

HTRF cAMP Gs HiRange kit - 100K pts - Lot n° : 21E

Accepted

Rejected

| | |
|-----------------------------------|----------------------------|
| Quality control date : 2024-10-22 | 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.