

HTRF® Research Reagents

HTRF cAMP Gi kit - 1K pts

Product Number : 62AM9PEB Lot Number : 06P

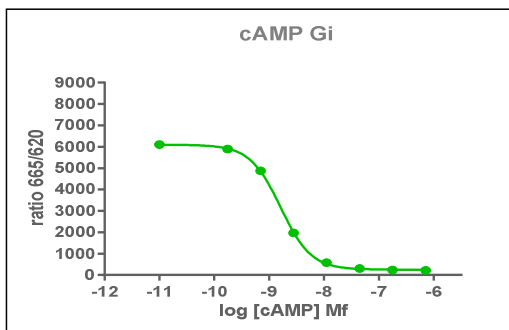
Expiration date : 2024-06-15 Document version : 1

| Reagents : | Lot n° |
|---|--------|
| HTRF Lysis & Detect. Buf. 8 - 13 mL | 06RA |
| HTRF Stimulation Buf. 1 (5X) - 8 mL | 24RA |
| HTRF cAMP Gs Dynamic & Gi kit - Std | 17AR |
| HTRF cAMP Gi kit - 1K pts Eu Cryptate reagent | 06RA |
| HTRF cAMP Gi kit - 1K pts d2 antibody | 06RA |
| HTRF cAMP Gi kit - modop | 5 |

Results :

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

| cAMP | Conc. (nM) | DF % |
|-------|------------|------|
| Std 0 | 0 | 2784 |
| Std 1 | 0.69 | 2683 |
| Std 2 | 2.78 | 2200 |
| Std 3 | 11.12 | 829 |
| Std 4 | 44.5 | 174 |
| Std 5 | 178 | 44 |
| Std 6 | 712 | 12 |
| Std 7 | 2848 | 1 |



Assay characteristics :

| | Found value | Specifications |
|----------------|-------------|----------------|
| S/B | 29 | 27-40 |
| Hillslope | -1.556 | -1.35-2.04 |
| IC50 Nm | 1.604 | 1.25-3 |
| blanc k 620 mm | 30156 | |

HTRF cAMP Gi kit - 1K pts - Lot n° : 06P

Accepted

Rejected

Quality control date : 2023-10-16 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.