

HTRF® Research Reagents

HTRF cGMP kit - 500 pts

Product Number : 62GMP2PEG Lot Number : 17D

Expiration date : 13/04/2025 Document version : 1

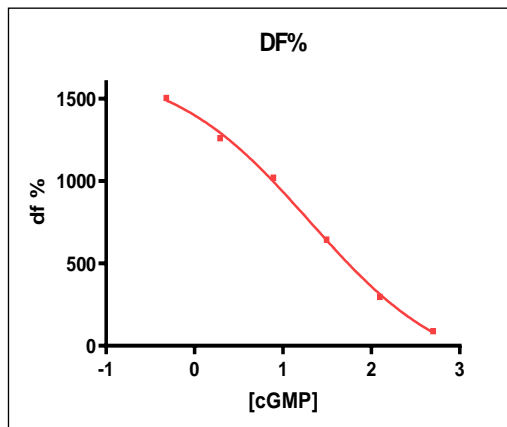
| Reagents : | Lot n° |
|---|--------|
| Diluent 1 - 20 mL | 19A |
| Lysis & Detection Buffer 1 - 7 mL | 14RA |
| cGMP kit - 500 tests d2 reagent | 17A |
| cGMP kit - modop | 6 |
| cGMP kit standard | 14B |
| cGMP kit - 500 tests Eu Cryptate antibody | 17A |

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)

Assay characteristics :

| cGMP | Conc. (nMf) | DF % |
|-------|-------------|-------|
| Std 0 | 0 | 1595% |
| Std 1 | 0,48 | 1504% |
| Std 2 | 1,95 | 1261% |
| Std 3 | 7,8 | 1020% |
| Std 4 | 31,5 | 644% |
| Std 5 | 120 | 296% |
| Std 6 | 500 | 89% |



Assay characteristics :

| | Found value | Specifications |
|------------------|-------------|----------------|
| DF max % | 1595% | 1450% - 2100% |
| EC 50 | 20 nMf | 10 nMf – 21nMf |
| Detection limite | 0.3 nMf | ≤ 1 nMf |

Accepted

Rejected

HTRF cGMP kit - 500 pts - Lot n° : 17D

Quality control date : 2023-09-15 Supervisor's initials : FMV

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.