

HTRF® Research Reagents

HTRF cGMP kit - 500 pts

Product Number : 62GMP2PEG Lot Number : 17F

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

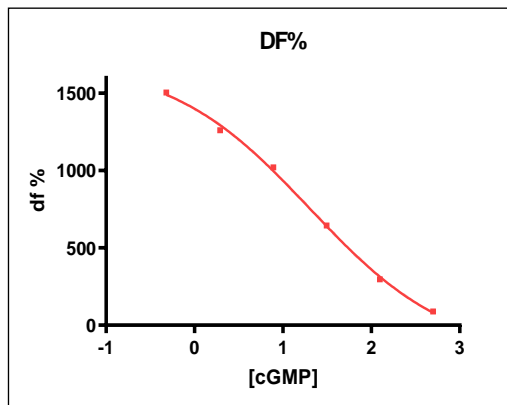
Reagents :	Lot n°
Diluent 1 - 20 mL	20RA
Lysis & Detection Buffer 1 - 7 mL	14RA
cGMP kit - 500 tests d2 reagent	17A
cGMP kit - modop	6
cGMP kit standard	14B
cGMP kit - 500 tests Eu Cryptate antibody	17A

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° : 17F

Quality control date : 2023-09-19 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.