

HTRF® Research Reagents

HTRF cGMP kit - 500 pts

Product Number : 62GM2PEG Lot Number : 20B

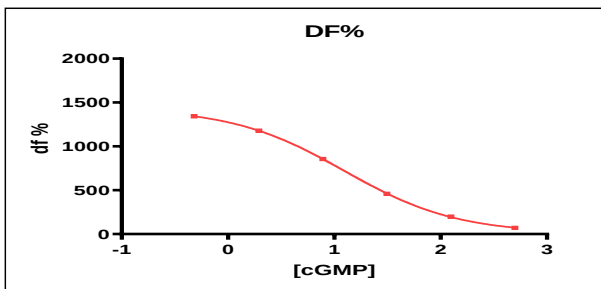
Expiration date : 2027-06-16 Document version : 1

Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20RA
HTRF Lysis & Detect. Buf. 1 - 7 mL	14RA
HTRF cGMP kit - 500 pts d2 reagent	20A
HTRF cGMP kit - Std	16A
HTRF cGMP kit - 500 pts Eu Cryptate antibody	20A

Results :

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

cGMP	Conc. (nMf)	DF %
Std 0	0	1504
Std 1	0,48	1342
Std 2	1,95	1179
Std 3	7,8	856
Std 4	31,5	459
Std 5	120	199
Std 6	500	70



Assay characteristics :

	Found value	Specifications Rev 01
DF max %	1504%	1450% - 2100%
EC 50	12,39 nMf	10 nMf - 21nMf
Detection limite	0,23nMf	≤ 1 nMf

HTRF cGMP kit - 500 pts - Lot n° : 20B

☒ Accepted

☐ Rejected

Quality control date : 2025-04-15 Supervisor's initials : PL

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.