

HTRF® Research Reagents

HTRF cGMP kit - 500 pts

Product Number : 62GM2PEG Lot Number : 19B

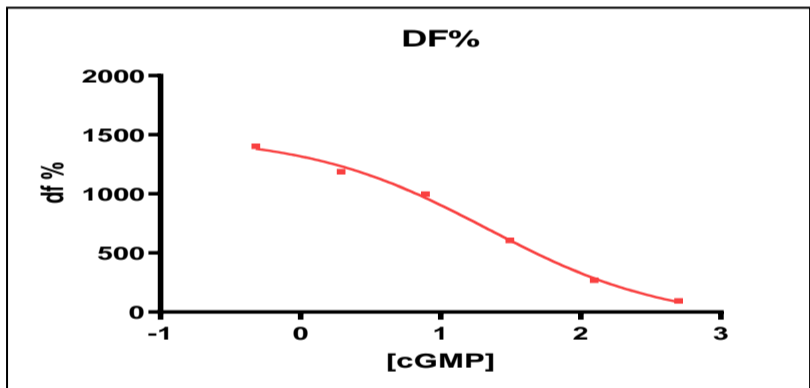
Expiration date : 2025-08-03 Document version : 1

Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20RA
HTRF Lysis & Detect. Buf. 1 - 7 mL	14RA
HTRF cGMP kit - 500 pts d2 reagent	19A
HTRF cGMP kit - Std	16A
HTRF cGMP kit - 500 pts Eu Cryptate antibody	17RA

Results :

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

cGMP	Conc. (nMf)	DF %
Std 0	0	1510%
Std 1	0,48	1403%
Std 2	1,95	1187%
Std 3	7,8	998%
Std 4	31,5	605%
Std 5	120	269%
Std 6	500	94%



Assay characteristics :

	Found value	Specifications
DF max %	1510%	1450% - 2100%
EC 50	21,5 nMf	10 nMf – 21nMf
Detection limite	0,55nMf	≤ 1 nMf

HTRF cGMP kit - 500 pts - Lot n° : 19C

Accepted

Rejected

Quality control date : 2025-01-27	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.