

HTRF® Research Reagents

HTRF cGMP kit - 10K pts

Product Number : 62GM2PEH Lot Number : 17G

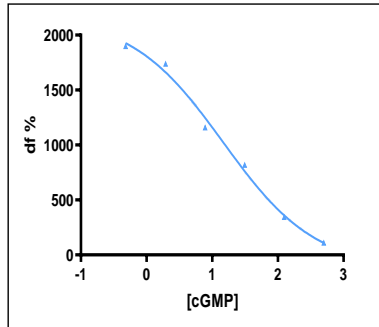
Expiration date : 2024-08-24 Document version : 1

Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20RA
HTRF Lysis & Detect. Buf. 2 - 105 mL	26RA
HTRF cGMP kit - 10K pts d2 reagent	17RB
HTRF cGMP kit - modop	6
HTRF cGMP kit - Std	14RB
HTRF cGMP kit - 10K pts Eu Cryptate antibody	17RA

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)

cGMP	Conc. (nMf)	DF %
Std 0	0	1942%
Std 1	0,488	1898%
Std 2	1,953	1738%
Std 3	7,81	1158%
Std 4	31,25	818%
Std 5	125	342%
Std 6	500	109%



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	0,89	<1 NmF
EC 50	14,7	10-21 NmF
DF% MAX	1942	1450-2100

HTRF cGMP kit - 10K pts - Lot n° : 17G

Accepted

Rejected

Quality control date : 2023-10-25 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.