

HTRF® Research Reagents

### HTRF cGMP kit - 500 pts

Product Number : 62GM2PEG Lot Number : 18B

Expiration date : 2025-06-16 Document version : 1

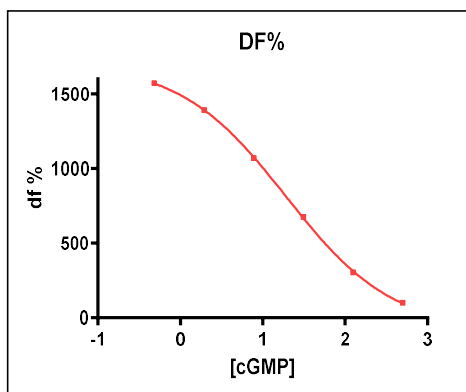
Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20RA
HTRF Lysis & Detect. Buf. 1 - 7 mL	14RA
HTRF cGMP kit - 500 pts d2 reagent	18AR
HTRF cGMP kit - modop	6
HTRF cGMP kit - Std	15B
HTRF cGMP kit - 500 pts Eu Cryptate antibody	17RA

**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)

**Assay characteristics :**

cGMP	Conc. (nMf)	DF %
Std 0	0	1704%
Std 1	0,48	1572%
Std 2	1,95	1392%
Std 3	7,8	1071%
Std 4	31,5	676%
Std 5	120	304%
Std 6	500	100%



HTRF cGMP kit - 500 pts - Lot n° : 18B

Accepted  Rejected

Quality control date : 2024-02-07	Supervisor's initials : AC
-----------------------------------	----------------------------

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.