

HTRF® Research Reagents

HTRF cAMP Gs HiRange kit - 1K pts

Product Number : 62AM6PEB Lot Number : 22B

Expiration date : 2026-04-14 Document version : 1

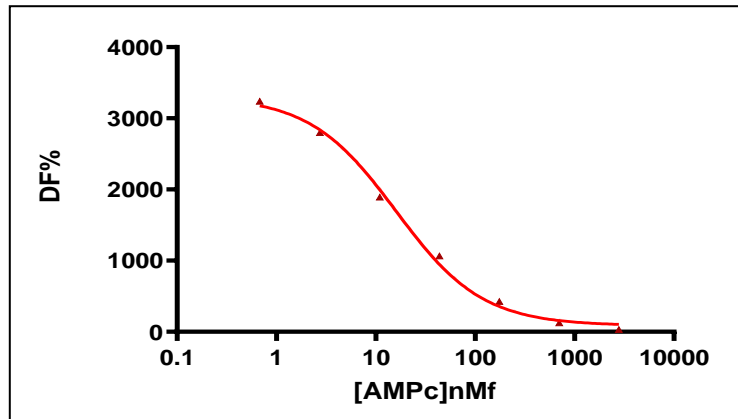
Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 1 - 13 mL	15A
HTRF Stimulation Buf. 1 (5X) - 8 mL	37A
HTRF cAMP Gs HiRange kit - Std	20A
HTRF cAMP Gs HiRange kit - 1K pts Eu	22A
HTRF cAMP Gs HiRange kit - 1K pts d2	22A

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 :

cAMP	Conc. (nMf)	DF (%)
Std 0	0	3626
Std 1	0,68	3372
Std 2	2,73	2856
Std 3	10,94	2030
Std 4	43,75	1040
Std 5	175	367
Std 6	700	116
Std 7	2800	22



Assay characteristics :

	Found value	Specifications Rev 01
DF % max	3656%	2900 % - 4150%
EC50	14,5	10 nMf - 22 nMf

HTRF cAMP Gs HiRange kit - 1K pts - Lot n° : 22B

Accepted

Rejected

Quality control date : 2025-01-20	Supervisor's initials : MP
-----------------------------------	----------------------------

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.