

HTRF® Research Reagents

HTRF cAMP Gs HiRange kit - 20K pts

Product Number : 62AM6PEC Lot Number : 24A

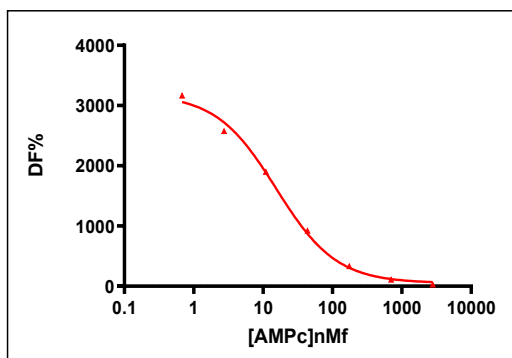
Expiration date : 2027-01-19 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 2 - 200 mL	33A
HTRF Stimulation Buf. 1 (5X) - 100 mL	46A
HTRF cAMP Gs HiRange kit - Std	21A
HTRF cAMP Gs HiRange kit - 20K pts Eu	23A
HTRF cAMP Gs HiRange kit - 20K pts d2	24A
/	/

Results :

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

cAMP	Conc. (nMf)	DF (%)
Std 0	0	3342
Std 1	0,68	3165
Std 2	2,73	2576
Std 3	10,94	1899
Std 4	43,75	923
Std 5	175	333
Std 6	700	109
Std 7	2800	28



Assay characteristics :

	Found value	Specifications Rev 01
DF% max	3342%	2900 % - 4150%
EC 50	15,43 nMf	10 nMf - 22 nMf

HTRF cAMP Gs HiRange kit - 20K pts - Lot n° : 24A

☒ Accepted☐ Rejected

Quality control date : 2025-08-27 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.