

HTRF® Research Reagents

HTRF cAMP Gs HiRange kit - 20K pts

Product Number : 62AM6PEC Lot Number : 22E

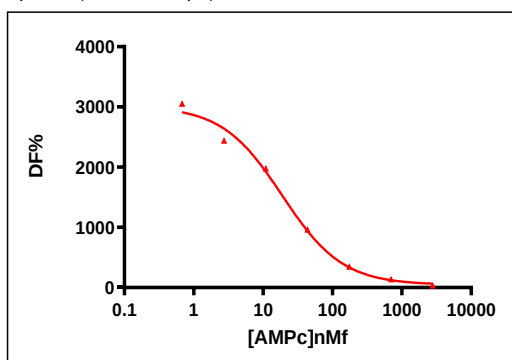
Expiration date : 2026-11-25 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 2 - 200 mL	32A
HTRF Stimulation Buf. 1 (5X) - 100 mL	44A
HTRF cAMP Gs HiRange kit - Std	20A
HTRF cAMP Gs HiRange kit - 20K pts Eu	22A
HTRF cAMP Gs HiRange kit - 20K pts d2	22A
/	/

Results :

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

cAMP	Conc. (nMf)	DF (%)
Std 0	0	3252
Std 1	0,68	3053
Std 2	2,73	2437
Std 3	10,94	1975
Std 4	43,75	963
Std 5	175	344
Std 6	700	134
Std 7	2800	30



Assay characteristics :

	Found value	Specifications Rev 01
DF% max	3252%	2900 % - 4150%
EC 50	16,3 nMf	10 nMf - 22 nMf

HTRF cAMP Gs HiRange kit - 20K pts - Lot n° : 22E

☒ Accepted☐ Rejected

Quality control date : 2025-06-18 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.