

HTRF® Research Reagents

HTRF cAMP Gs HiRange kit - 20K pts

Product Number : 62AM6PEC Lot Number : 21A

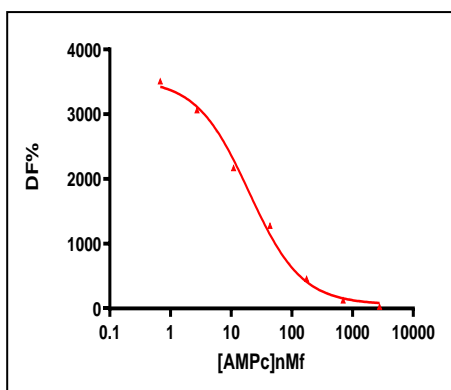
Expiration date : 2025-10-15 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 2 - 200 mL	30A
HTRF Stimulation Buf. 1 (5X) - 100 mL	31A
HTRF cAMP Gs HiRange kit - Std	19A
HTRF cAMP Gs HiRange kit - 20K pts Eu	21A
HTRF cAMP Gs HiRange kit - 20K pts d2	21A

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	3758
Std 1	0,68	3509
Std 2	2,73	3063
Std 3	10,94	2167
Std 4	43,75	1278
Std 5	175	458
Std 6	700	124
Std 7	2800	29



Assay characteristics :

	Found value	Specifications Rev 01
DF% max	3758%	2900 % - 4150%
EC 50	17,42	10 nMf - 22 nMf

HTRF cAMP Gs HiRange kit - 20K pts - Lot n° : 21A

Accepted

Rejected

Quality control date : 2024-05-17	Supervisor's initials : FL
-----------------------------------	----------------------------

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.