

HTRF® Research Reagents

HTRF cAMP Gs HiRange kit - 20K pts

Product Number : 62AM6PEC Lot Number : 20H

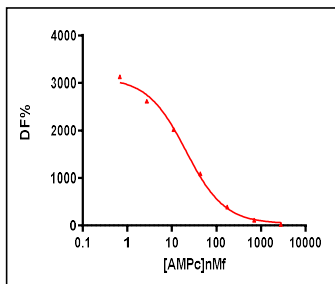
Expiration date : 2025-06-13 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 2 - 200 mL	28AR
HTRF Stimulation Buf. 1 (5X) - 100 mL	28AR
HTRF cAMP Gs HiRange kit - Std	16RA
HTRF cAMP Gs HiRange kit - 20K pts Eu	20RA
HTRF cAMP Gs HiRange kit - 20K pts d2	20RA
/	/

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	3284
Std 1	0,68	3127
Std 2	2,73	2616
Std 3	10,94	2019
Std 4	43,75	1088
Std 5	175	385
Std 6	700	112
Std 7	2800	25



Assay characteristics :

	Found value	Specifications Rev 01
DF% max	3284%	2900 % - 4150%
EC 50	19,19	10 nMf - 22 nMf
	/	/
	/	/

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

Accepted

Rejected

Quality control date : 2024-01-17 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.