

HTRF® Research Reagents

HTRF cAMP Gs Dynamic kit - 20K pts

Product Number : 62AM4PEC Lot Number : 35C

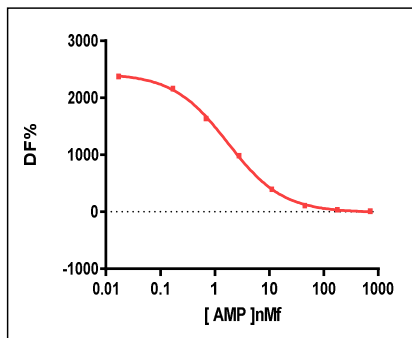
Expiration date : 2025-04-30 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 2 - 200 mL	28RA
HTRF Stimulation Buf. 1 (5X) - 100 mL	27AR
HTRF cAMP Gs Dynamic & Gi kit - Std	17AR
HTRF cAMP Gs Dynamic kit - 20K pts Eu	35AR
HTRF cAMP Gs Dynamic kit - 20K pts d2	34BR
HTRF cAMP Gs Dynamic kit - modop	8

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)

cAMP	Conc. (nMf)	DF %
Std 0	0	2375
Std 1	0.17	2162
Std 2	0.69	1635
Std 3	2.78	984
Std 4	11.12	390
Std 5	44.5	107
Std 6	178	34
Std 7	712	8



Assay characteristics :

	Found value	Specifications Rev 01
DF % max	2375%	1700 - 2700 %
S/B	22.8	>15
EC 50	1.7	< ou = 5.6 nMf
	/	/

HTRF cAMP Gs Dynamic kit - 20K pts - Lot n° : 35C

Accepted

Rejected

Quality control date : 2023-11-30	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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.