

HTRF® Research Reagents

HTRF cAMP Gs Dynamic kit - 20K pts

Product Number : 62AM4PEC Lot Number : 36B

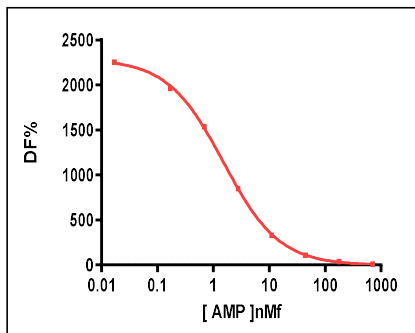
Expiration date : 2025-06-13 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 2 - 200 mL	28AR
HTRF Stimulation Buf. 1 (5X) - 100 mL	28AR
HTRF cAMP Gs Dynamic & Gi kit - Std	17AR
HTRF cAMP Gs Dynamic kit - 20K pts Eu	36AR
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	2256
Std 1	0,17	1964
Std 2	0,69	1536
Std 3	2,78	850
Std 4	11,12	328
Std 5	44,5	110
Std 6	178	37
Std 7	712	9



Assay characteristics :

	Found value	Specifications Rev 01
DF % max	2256%	1700 - 2700 %
S/B	21,5	>15
EC 50	1,5	< ou = 5.6 nMf
	/	/

HTRF cAMP Gs Dynamic kit - 20K pts - Lot n° : 36B

Accepted Rejected

Quality control date : 2024-01-15 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.