

HTRF® Research Reagents

HTRF cAMP Gs Dynamic kit - 20K pts

Product Number : 62AM4PEC Lot Number : 34J

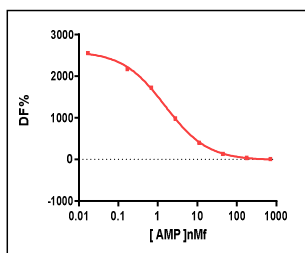
Expiration date : 2025-04-10 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 2 - 200 mL	28RA
HTRF Stimulation Buf. 1 (5X) - 100 mL	26RA
HTRF cAMP Gs Dynamic & Gi kit - Std	17AR
HTRF cAMP Gs Dynamic kit - 20K pts Eu	34BR
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	2557
Std 1	0.17	2175
Std 2	0.69	1729
Std 3	2.78	985
Std 4	11.12	406
Std 5	44.5	135
Std 6	178	39
Std 7	712	11



Assay characteristics :

	Found value	Specifications Rev 01
DF % max	2124%	1700 - 2700 %
S/B	20	>15
EC 50	1.9	< ou = 5.6 nMf
	/	/

HTRF cAMP Gs Dynamic kit - 20K pts - Lot n° : 34J

Accepted

Rejected

Quality control date : 2023-10-24 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.