

HTRF® Research Reagents

HTRF cAMP Gs Dynamic kit - 1K pts

Product Number : 62AM4PEB Lot Number : 16E

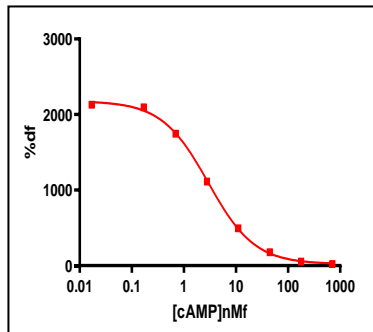
Expiration date : 2026-03-31 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 1 - 13 mL	14AR
HTRF Stimulation Buf. 1 (5X) - 8 mL	36A
HTRF cAMP Gs Dynamic & Gi kit - Std	18A
HTRF cAMP Gs Dynamic kit - 1K pts Eu	16A
HTRF cAMP Gs Dynamic kit - 1K pts d2	16A

Results :

Incubation : 1H A RT
 Plate : PROXIPLATE 384 WELLS WHITE PLATE SV
 Reader : Envision
 Fitting model : non linear regression, sigmoidal dose reponse (variable slope)

cAMP	Conc. (nMf)	DF %
Std 0	0	2131
Std 1	0,17	2096
Std 2	0,69	1746
Std 3	2,78	1113
Std 4	11,12	494
Std 5	44,5	177
Std 6	178	53
Std 7	712	20



Assay characteristics :

	Found value	Specifications Rev 01
DF MAX %	2131	1700-2700%
EC 50	2,9	< OU =5.6 nMf
	/	/
	/	/

HTRF cAMP Gs Dynamic kit - 1K pts - Lot n° : 16E

Accepted

Rejected

Quality control date : 2025-01-08	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.