

HTRF® Research Reagents

HTRF cAMP Gs Dynamic kit - 1K pts

Product Number : 62AM4PEB Lot Number : 16B

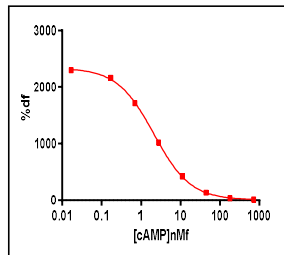
Expiration date : 10/15/2025 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 1 - 13 mL	14AR
HTRF Stimulation Buf. 1 (5X) - 8 mL	31A
HTRF cAMP Gs Dynamic & Gi kit - Std	17AR
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	2295
Std 1	0,17	2161
Std 2	0,69	1717
Std 3	2,78	1018
Std 4	11,12	423
Std 5	44,5	135
Std 6	178	39
Std 7	712	16



Assay characteristics :

	Found value	Specifications Rev 01
DF MAX %	2295	1700-2700%
EC 50	2,1	< OU =5.6 nMf
	/	/
	/	/

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

Accepted

Rejected

Quality control date : 2024-06-20	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.