

HTRF® Research Reagents

HTRF cAMP Gs Dynamic kit - 1K pts

Product Number : 62AM4PEB Lot Number : 15L

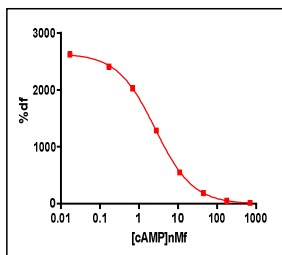
Expiration date : 2024-09-21 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 1 - 13 mL	14A
HTRF Stimulation Buf. 1 (5X) - 8 mL	24RA
HTRF cAMP Gs Dynamic & Gi kit - Std	17AR
HTRF cAMP Gs Dynamic kit - 1K pts Eu	15A
HTRF cAMP Gs Dynamic kit - 1K pts d2	15A
HTRF cAMP Gs Dynamic kit - modop	8

Results :

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

cAMP	Conc. (nMf)	DF %
Std 0	0	2626
Std 1	0.17	2412
Std 2	0.69	2031
Std 3	2.78	1289
Std 4	11.12	551
Std 5	44.5	190
Std 6	178	50
Std 7	712	17



Assay characteristics :

	Found value	Specifications Rev 01
DF MAX %	2626	1700-2700%
EC 50	2.5	< OU =5.6 nMf
S/B	/	/
	/	/

HTRF cAMP Gs Dynamic kit - 1K pts - Lot n° : 15L

Accepted

Rejected

Quality control date : 2023-09-14	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.