

HTRF® Research Reagents

HTRF cAMP Gs Dynamic kit - 1K pts

Product Number : 62AM4PEB Lot Number : 17B

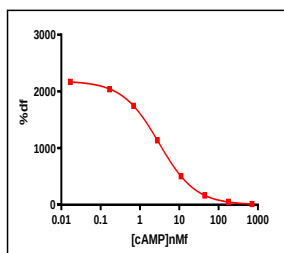
Expiration date : 2026-06-03 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 1 - 13 mL	15A
HTRF Stimulation Buf. 1 (5X) - 8 mL	39A
HTRF cAMP Gs Dynamic & Gi kit - Std	18A
HTRF cAMP Gs Dynamic kit - 1K pts Eu	17A
HTRF cAMP Gs Dynamic kit - 1K pts d2	17A
/	/
/	/

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	2169
Std 1	0,17	2038
Std 2	0,69	1746
Std 3	2,78	1138
Std 4	11,12	503
Std 5	44,5	164
Std 6	178	52
Std 7	712	13



Assay characteristics :

	Found value	Specifications Rev 01
DF MAX %	2169	1700-2700%
EC 50	3	< OU =5.6 nMf
/	/	/
/	/	/

HTRF cAMP Gs Dynamic kit - 1K pts - Lot n° : 17B

☒ Accepted☐ Rejected

Quality control date : 2025-03-19 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.