

HTRF® Research Reagents

HTRF cAMP Gi kit - 1K pts

Product Number : 62AM9PEB Lot Number : 09B

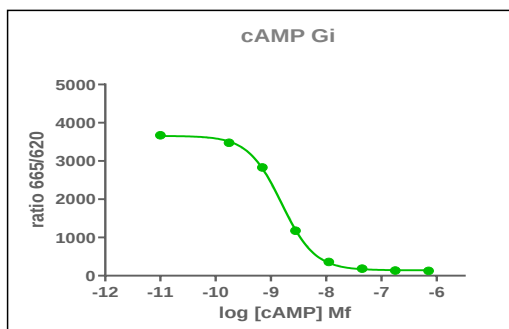
Expiration date : 2026-09-25 Document version : 1

Reagents :		Lot n°
HTRF Lysis & Detect. Buf. 8 - 13 mL		07A
HTRF Stimulation Buf. 1 (5X) - 8 mL		43A
HTRF cAMP Gs Dynamic & Gi kit - Std		19A
HTRF cAMP Gi kit - 1K pts Eu Cryptate reagent		09A
HTRF cAMP Gi kit - 1K pts d2 antibody		09A
/		/
/		/

Results :

Incubation : 1h at RT
Plate : 384 wells small volume white plate 20 µL
Reader : Envision
Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

cAMP	Conc. (nM)	DF %
Std 0	0	3178
Std 1	0,69	3006
Std 2	2,78	2429
Std 3	11,12	949
Std 4	44,5	223
Std 5	178	64
Std 6	712	20
Std 7	2848	12



Assay characteristics :

	Found value	Specifications
S/B	29	27-40
Hillslope	-1,44	-1,35-2,04
IC50 Nm	1,5	1.25-3
blanc k 620 mm	247600	

HTRF cAMP Gi kit - 1K pts - Lot n° : 09b

☒ Accepted☐ Rejected

Quality control date : 2025-06-17 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.