

HTRF® Research Reagents

HTRF cAMP Gi kit - 20K pts

Product Number : 62AM9PEC

Lot Number : 10A

Expiration date : 2027-01-19

Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 8 - 200 mL	07A
HTRF Stimulation Buf. 1 (5X) - 100 mL	46A
HTRF cAMP Gs Dynamic & Gi kit - Std	19A
HTRF cAMP Gi kit - 20K pts Eu Cryptate reagent	10A
HTRF cAMP Gi kit - 20K pts d2 antibody	10A
/	/
/	/

Results :

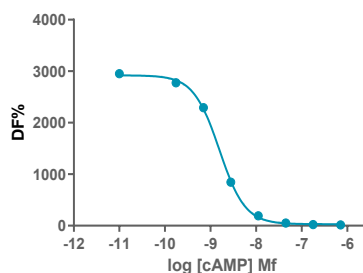
Incubation : 1h at RT

Plate : 384 wells small volume white plate 20 µL

Reader : Envision

Fitting model :

cAMP	Conc. (nMf)	DF %
Std 0	0	2950
Std 1	0,17	2775
Std 2	0,69	2293
Std 3	2,78	846
Std 4	11,12	192
Std 5	44,5	52
Std 6	178	22
Std 7	712	14



Assay characteristics :

	Found value	Specifications Rev 01
S/B	27	27-40
Hillslope	-1,54	between -1.35 and -2.04
IC50	1,55	between 1.25 and 3
blanc K 620 nm	161865	/

HTRF cAMP Gi kit - 20K pts - Lot n° : 10A

☒ Accepted☐ Rejected

Quality control date : 2025-09-25

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.