

HTRF® Research Reagents

HTRF cAMP Gi kit - 100K pts

Product Number : 62AM9PEJ Lot Number : 07B

Expiration date : 2025-04-30 Document version : 1

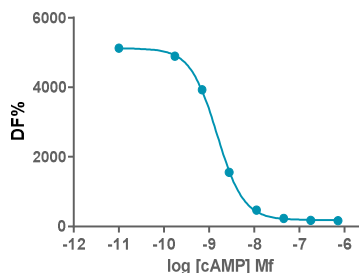
Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 8 - 200 mL	06RA
HTRF Stimulation Buf. 1 (5X) - 100 mL	27AR
HTRF cAMP Gs Dynamic & Gi kit - Std	17AR
HTRF cAMP Gi kit - 20K pts Eu Cryptate reagent	07RA
HTRF cAMP Gi kit - 20K pts d2 antibody	07RA
HTRF cAMP Gi kit - modop	5

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus

Fitting model :

cAMP	Conc. (nMf)	DF %
Std 0	0	3415
Std 1	0.17	3254
Std 2	0.69	2590
Std 3	2.78	968
Std 4	11.12	224
Std 5	44.5	58
Std 6	178	17
Std 7	712	9



Assay characteristics :

	Found value	Specifications Rev 01
S/B	32	27-40
Hillslope	-1.472	-1.35 ... -2.04
IC50	1.484	1.25-3
blanc K 620 nm	79216	/

HTRF cAMP Gi kit - 100K pts - Lot n° : 07B

Accepted

Rejected

Quality control date : 2023-11-30	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.