

HTRF® Research Reagents

HTRF cAMP Gi kit - 20K pts

Product Number : 62AM9PEC Lot Number : 06F

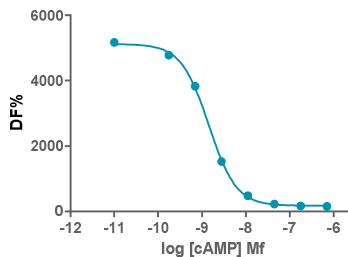
Expiration date : 2024-11-02 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 8 - 200 mL	06RA
HTRF Stimulation Buf. 1 (5X) - 100 mL	24RA
HTRF cAMP Gs Dynamic & Gi kit - Std	16RA
HTRF cAMP Gi kit - 20K pts Eu Cryptate reagent	06A
HTRF cAMP Gi kit - 20K pts d2 antibody	06A
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	3681
Std 1	0.17	3395
Std 2	0.69	2702
Std 3	2.78	1017
Std 4	11.12	251
Std 5	44.5	65
Std 6	178	21
Std 7	712	11



Assay characteristics :

	Found value	Specifications Rev 01
S/B	34	27-40
Hillslope	-1.385	between -1.35 and -2.04
IC50	1.418	between 1.25 and 3
blanc K 620 nm	94136	/

HTRF cAMP Gi kit - 20K pts - Lot n° : 06F

Accepted

Rejected

Quality control date : 2023-09-15	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.