

HTRF® Research Reagents

HTRF cAMP Gs HiRange kit - 100K pts

Product Number : 62AM6PEJ Lot Number : 20G

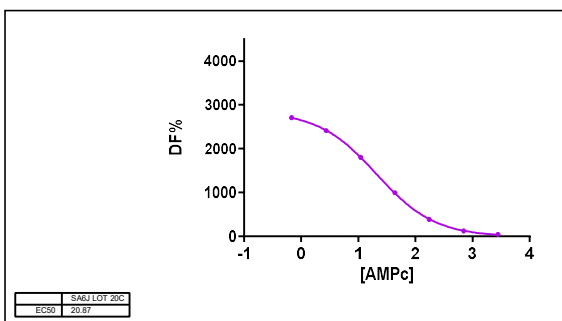
Expiration date : 2025-07-16 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 2 - 200 mL	28AR
HTRF Stimulation Buf. 1 (5X) - 100 mL	29AR
HTRF cAMP Gs HiRange kit - Std	19A
HTRF cAMP Gs HiRange kit - 20K pts Eu	20RA
HTRF cAMP Gs HiRange kit - 20K pts d2	20RA
/	/

Results :

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

cAMP	Conc. (nMf)	DF (%)
Std 0	0	2910
Std 1	0,68	2704
Std 2	2,73	2410
Std 3	10,94	1802
Std 4	43,75	990
Std 5	175	385
Std 6	700	123
Std 7	2800	42



Assay characteristics :

	Found value	Specifications Rev 01
DF% max	2910%	2900 % - 4150%
EC 50	20,8	10 nMf - 22 nMf
	/	/
	/	/

HTRF cAMP Gs HiRange kit - 100K pts - Lot n° : 20G

Accepted

Rejected

Quality control date : 2024-02-16	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.