

HTRF® Research Reagents

### HTRF cAMP Gs HiRange kit - 1K pts

Product Number : 62AM6PEB Lot Number : 15H

Expiration date : 2025-04-10 Document version : 1

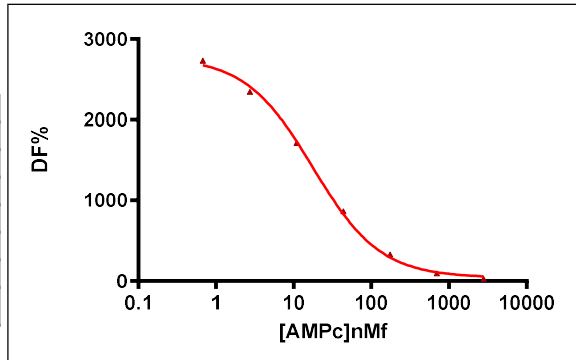
Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 1 - 13 mL	14AR
HTRF Stimulation Buf. 1 (5X) - 8 mL	26AR
HTRF cAMP Gs HiRange kit - Std	19A
HTRF cAMP Gs HiRange kit - 1K pts Eu	15RA
HTRF cAMP Gs HiRange kit - 1K pts d2	15RA
/	/
/	/

**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)

**Assay characteristics :**

cAMP	Conc. (nMf)	DF (%)
Std 0	0	3012
Std 1	0,68	2732
Std 2	2,73	2343
Std 3	10,94	1711
Std 4	43,75	864
Std 5	175	328
Std 6	700	97
Std 7	2800	25



**Assay characteristics :**

	Found value	Specifications Rev 01
DF % max	3012%	2900 % - 4150%
EC50	16,3	10 nMf - 22 nMf
	/	/
	/	/

HTRF cAMP Gs HiRange kit - 1K pts - Lot n° : 15H

Accepted

Rejected

Quality control date : 2024-02-08 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.