

HTRF® Research Reagents

### HTRF cAMP Gs Dynamic kit - 20K pts

Product Number : 62AM4PEC Lot Number : 39B

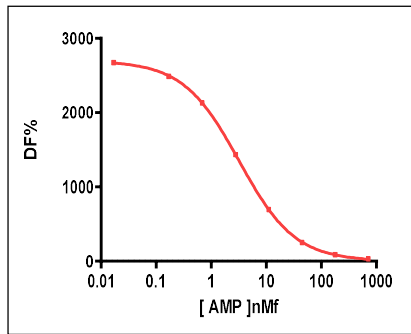
Expiration date : 2026-03-14 Document version : 1

Reagents :		Lot n°
HTRF Lysis & Detect. Buf. 2 - 200 mL		30A
HTRF Stimulation Buf. 1 (5X) - 100 mL		35A
HTRF cAMP Gs Dynamic & Gi kit - Std		17AR
HTRF cAMP Gs Dynamic kit - 20K pts Eu		39A
HTRF cAMP Gs Dynamic kit - 20K pts d2		39A

**Results :**

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

cAMP	Conc. (nMf)	DF %
Std 0	0	2673
Std 1	0,17	2485
Std 2	0,69	2132
Std 3	2,78	1433
Std 4	11,12	697
Std 5	44,5	250
Std 6	178	85
Std 7	712	30



**Assay characteristics :**

	Found value	Specifications Rev 01
DF % max	2673%	1700 - 2700 %
S/B	21,3	>15
EC 50	3,1	< ou = 5.6 nMf

HTRF cAMP Gs Dynamic kit - 20K pts - Lot n° : 39B

Accepted

Rejected

Quality control date : 2024-10-26	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.