

HTRF® Research Reagents

HTRF cAMP Gs Dynamic kit - 20K pts

Product Number : 62AM4PEC Lot Number : 40C

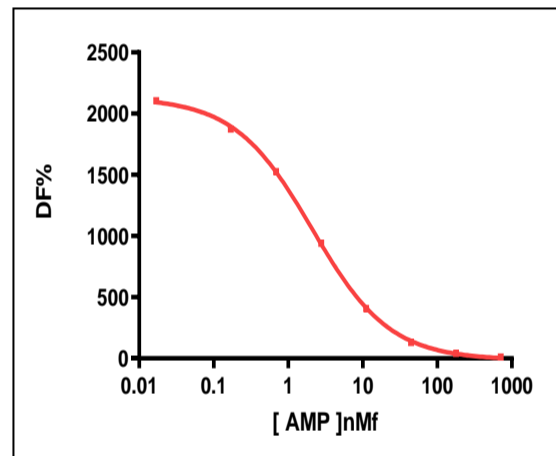
Expiration date : 2026-06-03 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 2 - 200 mL	31A
HTRF Stimulation Buf. 1 (5X) - 100 mL	39A
HTRF cAMP Gs Dynamic & Gi kit - Std	18A
HTRF cAMP Gs Dynamic kit - 20K pts Eu	40A
HTRF cAMP Gs Dynamic kit - 20K pts d2	40A

Results :

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

cAMP	Conc. (nMf)	DF %
Std 0	0	2105
Std 1	0,17	1873
Std 2	0,69	1527
Std 3	2,78	941
Std 4	11,12	407
Std 5	44,5	131
Std 6	178	41
Std 7	712	11



Assay characteristics :

	Found value	Specifications Rev 01
DF % max	2105%	1700 - 2700 %
S/B	19,9	>15
EC 50	2,1	< ou = 5.6 nMf

HTRF cAMP Gs Dynamic kit - 20K pts - Lot n° : 40C

Accepted

Rejected

Quality control date : 2025-02-19	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.