

HTRF® Research Reagents

HTRF cAMP Gs Dynamic kit - 1K pts

Product Number : 62AM4PEB Lot Number : 17E

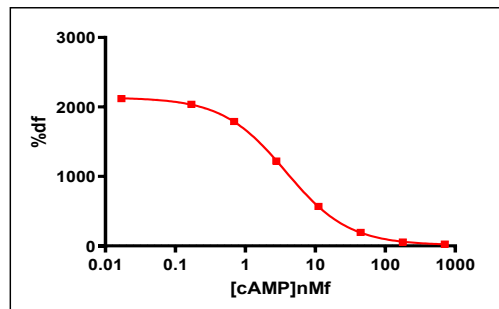
Expiration date : 2025-11-25 Document version :

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 1 - 13 mL	15A
HTRF Stimulation Buf. 1 (5X) - 8 mL	44A
HTRF cAMP Gs Dynamic & Gi kit - Std	19A
HTRF cAMP Gs Dynamic kit - 1K pts Eu	17A
HTRF cAMP Gs Dynamic kit - 1K pts d2	17A
/	/

Results :

Incubation : 1H A RT
Plate : PROXIPLATE 384 WELLS WHITE PLATE SV
Reader : Envision
Fitting model : non linear regression, sigmoidal dose response (variable slope)

cAMP	Conc. (nMf)	DF %
Std 0	0	2119
Std 1	0,17	2038
Std 2	0,69	1790
Std 3	2,78	1221
Std 4	11,12	568
Std 5	44,5	194
Std 6	178	56
Std 7	712	25



Assay characteristics :

	Found value	Specifications Rev 01
DF MAX %	2119	1700-2700%
EC 50	3,7	< OU =5.6 nMf

HTRF cAMP Gs Dynamic kit - 1K pts - Lot n° : 17E

☒ Accepted☐ Rejected

Quality control date : 2025-08-12	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.