

HTRF® Research Reagents

HTRF IP-One Gq kit - 1K pts

Product Number : 62IPAPEB Lot Number : 17D

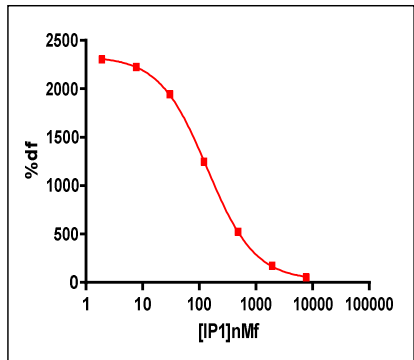
Expiration date : 2026-01-04 Document version : 1

Reagents :		Lot n°
HTRF Lysis & Detect. Buf. 5 - 13 mL		07RA
HTRF IP-One Gq kit - 1K pts Tb		17A
HTRF IP-One Gq kit - 1K pts d2		17A
HTRF IP-One Gq kit - modop		8
HTRF IP-One Gq kit - Std		15A
HTRF Stimulation Buf. 2 (5X) - 8 mL		37RA

Results :

Incubation : 1h at RT
 Plate : 384 wells regular white plate 30 µl
 Reader : Pherastar Plus
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

IP-One	Conc. (nMf)	DF %
Std 0	0	2294%
Std 1	1.9	2304%
Std 2	7.7	2225%
Std 3	30.1	1943%
Std 4	120.4	1248%
Std 5	481.6	525%
Std 6	1925	172%
Std 7	7700	51%



Assay characteristics :

	Found value	Specifications Rev 01
DF max %	2294%	1800 – 2600 %
EC 50	137.1 nMf	100 - 140 nMf
S/B	/	/
	/	/

HTRF IP-One Gq kit - 1K pts - Lot n° : 17D

Accepted

Rejected

Quality control date : 2023-10-02	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.