

HTRF® Research Reagents

HTRF IP-One Gq kit - 20K pts

Product Number : 62IPAPEC Lot Number : 22L

Expiration date : 2025-05-17 Document version : 1

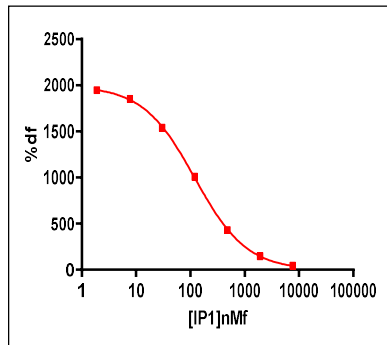
Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 6 - 200 mL	09RA
HTRF IP-One Gq kit - 20K pts Tb	22RB
HTRF IP-One Gq kit - 20K pts d2	22RB
HTRF IP-One Gq kit - modop	8
HTRF IP-One Gq kit - Std	16AR
HTRF Stimulation Buf. 2 (5X) - 100 mL	38AR
/	/

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)

Assay characteristics :

IP-One	Conc. (nMf)	DF %
Std 0	0	2018%
Std 1	1,9	1946%
Std 2	7,7	1852%
Std 3	30,1	1538%
Std 4	120,4	1008%
Std 5	481,6	430%
Std 6	1925	146%
Std 7	7700	43%



	Found value	Specifications Rev 01
DF MAX%	2018	1800-2600%
EC50	120,5	100-140 nMf
	/	/
	/	/

HTRF IP-One Gq kit - 20K pts - Lot n° : 22L  Accepted  Rejected

Quality control date : 2023-12-27 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.