

HTRF® Research Reagents

HTRF IP-One Gq kit - 20K pts

Product Number : 62IPAPEC Lot Number : 23B

Expiration date : 2026-08-02 Document version : 1

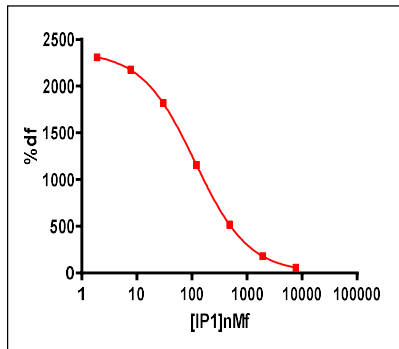
Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 6 - 200 mL	09RA
HTRF IP-One Gq kit - 20K pts Tb	23A
HTRF IP-One Gq kit - 20K pts d2	23A
HTRF IP-One Gq kit - Std	16AR
HTRF Stimulation Buf. 2 (5X) - 100 mL	38AR

Results :

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

Assay characteristics :

IP-One	Conc. (nMf)	DF %
Std 0	0	2325%
Std 1	1,9	2309%
Std 2	7,7	2175%
Std 3	30,1	1819%
Std 4	120,4	1156%
Std 5	481,6	517%
Std 6	1925	181%
Std 7	7700	55%



	Found value	Specifications Rev 01
DF MAX%	2325	1800-2600%
EC50	114,1	100-140 nMf
	/	/
	/	/

HTRF IP-One Gq kit - 20K pts - Lot n° : 23B

Accepted

Rejected

Quality control date : 2024-03-05 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.