

HTRF® Research Reagents

HTRF IP-One Gq kit - 100K pts

Product Number : 62IPAPEJ Lot Number : 22E

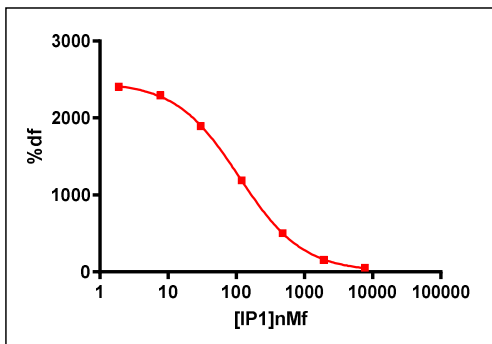
Expiration date : 2025-04-20 Document version : 1

Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 6 - 200 mL	09RA
HTRF IP-One Gq kit - 20K pts Tb	22RA
HTRF IP-One Gq kit - 20K pts d2	22RB
HTRF IP-One Gq kit - modop	8
HTRF IP-One Gq kit - Std	15RA
HTRF Stimulation Buf. 2 (5X) - 100 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	2399%
Std 1	1.9	2406%
Std 2	7.7	2293%
Std 3	30.1	1894%
Std 4	120.4	1189%
Std 5	481.6	502%
Std 6	1925	154%
Std 7	7700	50%



Assay characteristics :

	Found value	Specifications Rev 01
DF MAX%	2399	1800-2600%
EC50	110.8	100-140 nMf
	/	/
	/	/

HTRF IP-One Gq kit - 100K pts - Lot n° : 22E

Accepted

Rejected

Quality control date : 2023-10-16	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.