

HTRF® Research Reagents

HTRF IP-One Gq kit - 20K pts

Product Number : 62IPAPEC Lot Number : 24E

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

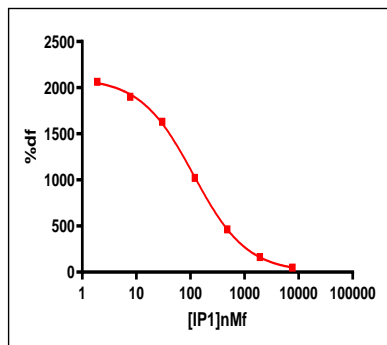
Reagents :	Lot n°
HTRF Lysis & Detect. Buf. 6 - 200 mL	10A
HTRF IP-One Gq kit - 20K pts Tb	24A
HTRF IP-One Gq kit - 20K pts d2	23A
HTRF IP-One Gq kit - Std	16AR
HTRF Stimulation Buf. 2 (5X) - 100 mL	40A
/	/
/	/

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	2071%
Std 1	1,9	2063%
Std 2	7,7	1901%
Std 3	30,1	1630%
Std 4	120,4	1022%
Std 5	481,6	464%
Std 6	1925	164%
Std 7	7700	48%



	Found value	Specifications Rev 01
DF MAX%	2071	1800-2600%
EC50	114,9	100-140 nMf
	/	/
	/	/

HTRF IP-One Gq kit - 20K pts - Lot n° : 24E Accepted Rejected

Quality control date : 2025-03-12 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.