

HTRF® Research Reagents

HTRF GTP Gi Bind. Kit - 100 pts

Product Number : 62GTPPET Lot Number : 05A

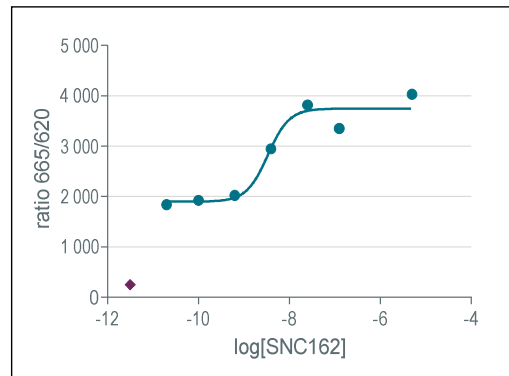
Expiration date : 2024-05-26 Document version : 1

Reagents :	Lot n°
HTRF Stimulation Buf. 3 - 10 mL	07AR
HTRF GTP Gi Bind. Kit - GDP	04AR
HTRF GTP Gi Bind. Kit - GTP gamma S	03RA
HTRF GTP Gi Bind. Kit - MgCL2	04AR
HTRF GTP Gi Bind. Kit - Gi protein control	02RA
HTRF GTP Gi Bind. Kit - 100 pts d2 antibody	05AR
HTRF GTP Gi Bind. Kit - 100 ptsEu Cryptate reagent	05AR
HTRF GTP Gi Bind. Kit - modop	2

Results :

Incubation : Overnight at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Pherastar Plus  
 Fitting model : log(agonist)vs.response. Variable slope (four parameters)

Agonist	Conc. (Mf)	ratio
Std 0	basal	1840
agonist 1	1.1E-10	1928
agonist 2	6.4E-10	2022
agonist 3	3.9E-09	2947
agonist 4	2.3E-08	3819
agonist 5	1.4E-07	3351
agonist 6	8.3E-07	3017
agonist 7	5.0E-06	4033
GTPyS		251
Gi protein		3792



Assay characteristics :

	Found value	Specifications Rev 01
S/B basal	7.3	3 - 8.5
S/B Gi control	17.3	15 - 25
EC50	3.4	3.3 - 10 nM
	/	/

HTRF GTP Gi Bind. Kit - 100 pts - Lot n° : 05A

Accepted

Rejected

Quality control date : 2023-08-18 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.