

HTRF® Research Reagents

HTRF GTP Gi Bind. Kit - 100 pts

Product Number : 62GTPPET Lot Number : 01D

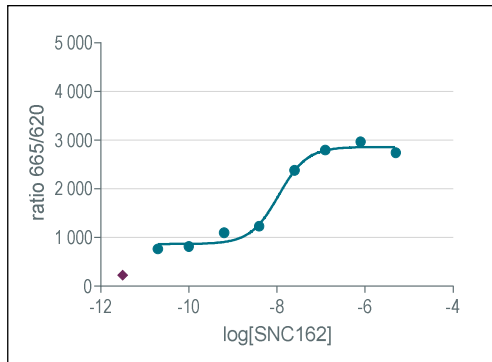
Expiration date : 2024-05-26 Document version : 1

Reagents :	Lot n°
HTRF Stimulation Buf. 3 - 10 mL	05RB
HTRF GTP Gi Bind. Kit - GDP	04AR
HTRF GTP Gi Bind. Kit - GTP gamma S	03RA
HTRF GTP Gi Bind. Kit - MgCL2	03AR
HTRF GTP Gi Bind. Kit - Gi protein control	02RA
HTRF GTP Gi Bind. Kit - 100 pts d2 antibody	01A
HTRF GTP Gi Bind. Kit - 100 ptsEu Cryptate reagent	01A
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	766
agonist 1	1.1E-10	814
agonist 2	6.4E-10	1096
agonist 3	3.9E-09	1232
agonist 4	2.3E-08	2383
agonist 5	1.4E-07	2796
agonist 6	8.3E-07	2967
agonist 7	5.0E-06	2742
GTPyS		222
Gi protein		3040



Assay characteristics :

	Found value	Specifications Rev 01
S/B basal	3.6	3 - 4.4
S/B Gi control	17.3	15 - 17.4
EC50	10.9 nM	9 -25 nM
	/	/

HTRF GTP Gi Bind. Kit - 100 pts - Lot n° : 01D

Accepted

Rejected

Quality control date :	2023-09-19	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.