

HTRF® Research Reagents

HTRF GTP Gi Bind. Kit - 20K pts

Product Number : 62GTPPEC Lot Number : 05CC

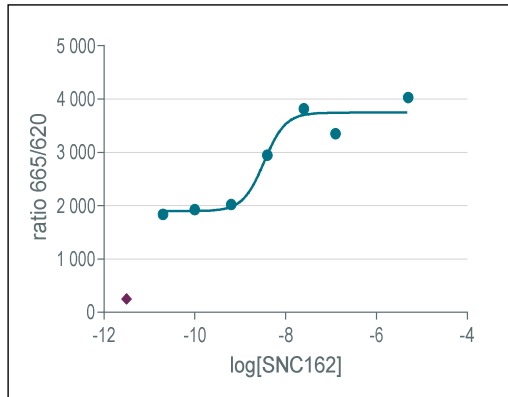
Expiration date : 2024-05-26 Document version : 1

Reagents :	Lot n°
HTRF GTP Gi Bind. Kit - 20K pts GDP	02RA
HTRF GTP Gi Bind. Kit - 20K pts d2 antibody	05AR
HTRF GTP Gi Bind. Kit - 20K pts GTP gamma S	04AR
HTRF GTP Gi Bind. Kit - 20K pts Eu Cryptate reagent	05AR
HTRF GTP Gi Bind. Kit - 20K pts MgCL2	04AR
HTRF GTP Gi Bind. Kit - Gi protein control	02RA
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 - 20K pts - Lot n° : 05CC

Accepted

Rejected

Quality control date : 2023-09-26	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.