

HTRF® Research Reagents

HTRF GTP Gi Bind. Kit - 500 pts

Product Number : 62GTPPEG Lot Number : 06D

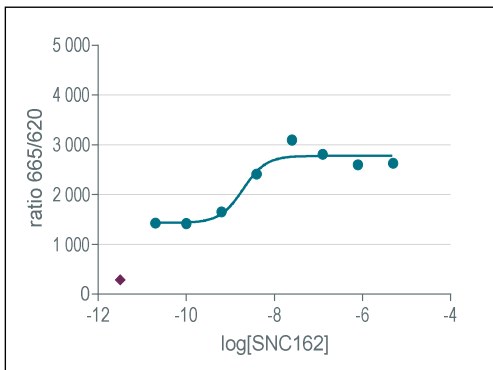
Expiration date : 2026-08-28 Document version : 1

Reagents :	Lot n°
HTRF Stimulation Buf. 3 - 10 mL	07AR
HTRF GTP Gi Bind. Kit - GDP	04AR
HTRF GTP Gi Bind. Kit - 500 pts d2 antibody	06AR
HTRF GTP Gi Bind. Kit - GTP gamma S	05AR
HTRF GTP Gi Bind. Kit - 500 pts Eu Cryptate reagent	06AR
HTRF GTP Gi Bind. Kit - MgCL2	04AR
HTRF GTP Gi Bind. Kit - Gi protein control	03AR
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	1429
agonist 1	1.1E-10	1418
agonist 2	6.4E-10	1657
agonist 3	3.9E-09	2413
agonist 4	2.3E-08	3097
agonist 5	1.4E-07	2809
agonist 6	8.3E-07	2604
agonist 7	5.0E-06	2633
/	/	/
GTPyS	/	291
Gi protein	/	3737



Assay characteristics :

	Found value	Specifications Rev 01
S/B basal	4.9	3 - 8.5
S/B Gi control	18.5	15 - 25
EC50	1.9	3.3 -10 nM
	/	/

Batch validated with EC50 a little bit lower than specification

HTRF GTP Gi Bind. Kit - 500 pts - Lot n° : 06D

Accepted

Rejected

Quality control date :	2023-10-23	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.