

HTRF® Research Reagents

HTRF GTP Gi Bind. Kit - 500 pts

Product Number : 62GTPPEG

Lot Number : 09J

Expiration date : 2026-11-08

Document version : 1

Reagents :	Lot n°
HTRF Stimulation Buf. 3 - 10 mL	12AR
HTRF GTP Gi Bind. Kit - GDP	10A
HTRF GTP Gi Bind. Kit - 500 pts d2 antibody	09AR
HTRF GTP Gi Bind. Kit - GTP gamma S	08A
HTRF GTP Gi Bind. Kit - 500 pts Eu Cryptate reagent	09AR
HTRF GTP Gi Bind. Kit - MgCL2	05A
HTRF GTP Gi Bind. Kit - Gi protein control	04AR
/	/
/	/

Results :

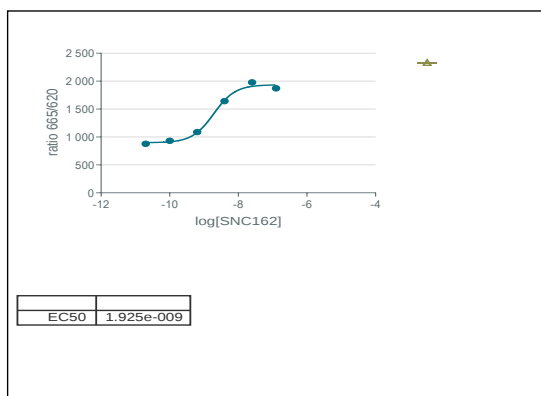
Incubation : Overnight at RT

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	879
agonist 1	1,1E-10	932
agonist 2	6,4E-10	1088
agonist 3	3,9E-09	1642
agonist 4	2,3E-08	1979
agonist 5	1,4E-07	1872
agonist 6	8,3E-07	1811
agonist 7	5,0E-06	1773
/	/	/
GTPyS	/	169
Gi protein	/	2906



Assay characteristics :

	Found value	Specifications Rev 01
S/B basal	5,2	3 - 8.5
S/B Gi control	18,9	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° : 09J

☒ Accepted☐ Rejected

Quality control date : 2025-04-04

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.