

HTRF® Research Reagents

HTRF GTP Gi Bind. Kit - 20K pts

Product Number : 62GTPPEC Lot Number : 09F

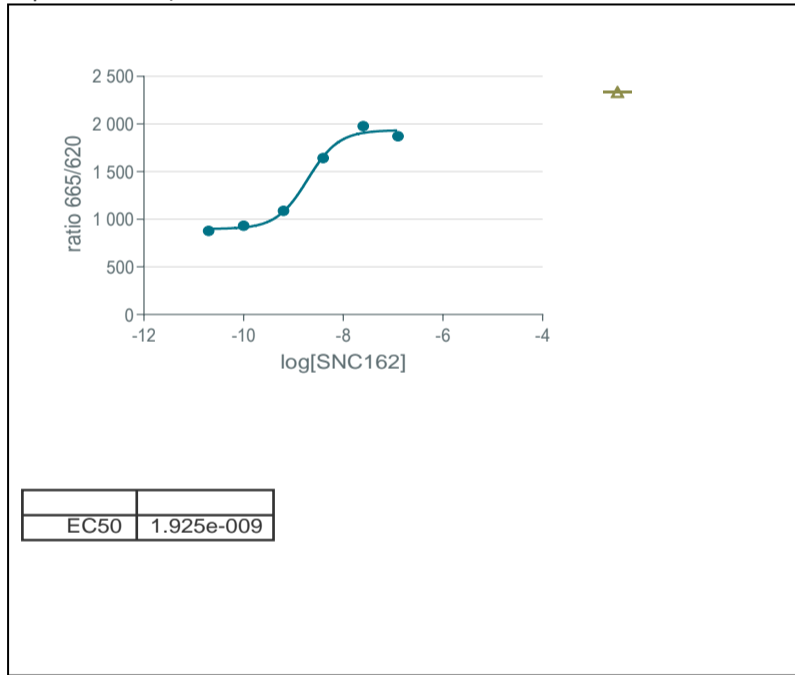
Expiration date : 2026-10-18 Document version : 1

Reagents :	Lot n°
HTRF Stimulation Buf. 3 - 200 mL	13A
HTRF GTP Gi Bind. Kit - 20K pts GDP	06AR
HTRF GTP Gi Bind. Kit - 20K pts d2 antibody	09AR
HTRF GTP Gi Bind. Kit - 20K pts GTP gamma S	08A
HTRF GTP Gi Bind. Kit - 20K pts Eu Cryptate reagent	09AR
HTRF GTP Gi Bind. Kit - 20K pts MgCL2	05AR
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	/	172
Gi protein	/	2793



Assay characteristics :

	Found value	Specifications Rev 01
S/B basal	5,1	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 - 20K pts - Lot n° : 09F

Accepted

Rejected

Quality control date : 2024-12-16 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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.