

HTRF® Research Reagents

HTRF GTP Gi Bind. Kit - 500 pts

Product Number : 62GTPPEG Lot Number : 09A

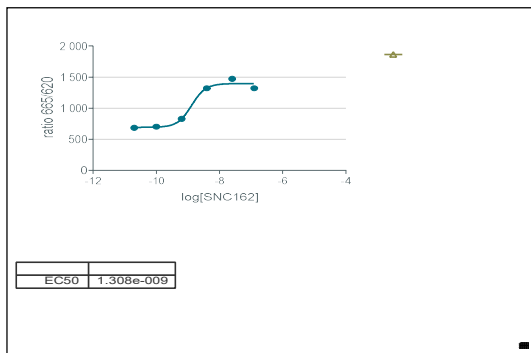
Expiration date : 2026-08-28 Document version : 1

Reagents :	Lot n°
HTRF Stimulation Buf. 3 - 10 mL	11AR
HTRF GTP Gi Bind. Kit - GDP	05AR
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	09AR
HTRF GTP Gi Bind. Kit - MgCL2	04AR
HTRF GTP Gi Bind. Kit - Gi protein control	04AR
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	686
agonist 1	1,1E-10	680
agonist 2	6,4E-10	831
agonist 3	3,9E-09	1320
agonist 4	2,3E-08	1474
agonist 5	1,4E-07	1322
agonist 6	8,3E-07	1344
agonist 7	5,0E-06	1303
/	/	/
GTPyS	/	171
Gi protein	/	2793



Assay characteristics :

	Found value	Specifications Rev 01
S/B basal	4	3 - 8.5
S/B Gi control	18,9	15 - 25
EC50	1,3	3.3 - 10 nM
	/	/

Batch validated with EC50 a little bit lower than specification

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

Accepted

Rejected

Quality control date : 2023-11-17 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.