

HTRF® Research Reagents

HTRF GTP Gi Bind. Kit - 100 pts

Product Number : 62GTPPET Lot Number : 05E

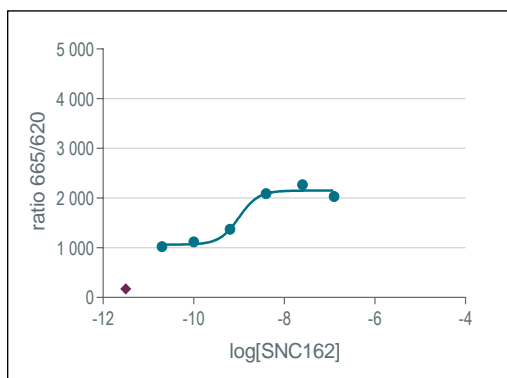
Expiration date : 2026-07-26 Document version : 1

Reagents :	Lot n°
HTRF Stimulation Buf. 3 - 10 mL	12AR
HTRF GTP Gi Bind. Kit - GDP	10A
HTRF GTP Gi Bind. Kit - GTP gamma S	07A
HTRF GTP Gi Bind. Kit - MgCL2	07A
HTRF GTP Gi Bind. Kit - Gi protein control	04AR
HTRF GTP Gi Bind. Kit - 100 pts d2 antibody	05AR
HTRF GTP Gi Bind. Kit - 100 ptsEu Cryptate reagent	05AR
/	/
/	/

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	1022
agonist 1	1,1E-10	1118
agonist 2	6,4E-10	1371
agonist 3	3,9E-09	2090
agonist 4	2,3E-08	2268
agonist 5	1,4E-07	2033
agonist 6	8,3E-07	1944
agonist 7	5,0E-06	1944
GTPyS		174
Gi protein		3576



Assay characteristics :

	Found value	Specifications Rev 01
S/B basal	5,9	3 - 8.5
S/B Gi control	24,1	15 - 25
EC50	9,9	3.3 -10 nM
/	/	/

HTRF GTP Gi Bind. Kit - 100 pts - Lot n° : 05E

Accepted

Rejected

Quality control date : 2025-03-11 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.