

HTRF® Research Reagents

### HTRF GST check kit - 500 pts

Product Number : 62GSTPEG Lot Number : 11C

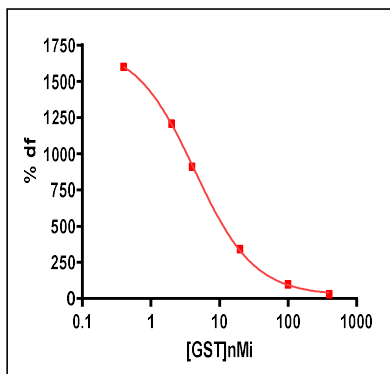
Expiration date : 2025-06-16 Document version : 1

Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20RA
HTRF Reconstitution Buf. 2 - 13 mL	13RA
HTRF GST check kit - 500 pts Eu Cryptate antibody	11RA
HTRF GST check kit - 500 pts XL665 reagent	10RA
HTRF GST check kit - Std	09RA

**Results :**

Incubation : 2h at RT  
 Plate : 384 wells small volume white plate 20 µL (Greiner)  
 Reader : Pherastar Plus  
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

GST	Conc. (nMi)	DF %
Std 0	0	1717%
Std 1	0,4	1601%
Std 2	2	1208%
Std 3	4	910%
Std 4	20	342%
Std 5	100	98%
Std 6	400	30%



**Assay characteristics :**

	Found value	Specifications Rev 03
Detection limit	0,22	< 0.26 nM
EC 50	4,31	4-10 nM
DF max	1717	1350-1800%

HTRF GST check kit - 500 pts - Lot n° : 11C

Accepted

Rejected

Quality control date : 2024-10-02	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.