

HTRF® Research Reagents

HTRF STING p-S365 (m) kit - 500 pts

Product Number : 64MSTGPEG Lot Number : 01J

Expiration date : 2025-03-08 Document version : 1

Reagents :	Lot n°
HTRF STING p-S365 (m) kit - 500 pts d2 antibody	01RA
HTRF STING p-S365 (m) kit - 500 pts Eu Cryptate antibody	01RA
HTRF STING p-S365 (m) kit - Ctrl lysate	02RA
HTRF STING p-S365 (m) kit - modop	1
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	14A
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28A
HTRF P-T prot. - Detect. Buf. 2 mL	32A

Results :

Incubation : 2h at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Pherastar Plus  
 Fitting model : /

Assay characteristics :

	Found value	Specifications Rev 01
Negative	407	/
Control Lysate	3858	/
S/B	9.5	>5
	/	/

HTRF STING p-S365 (m) kit - 500 pts - Lot n° : 01J

Accepted

Rejected

Quality control date : 2022-02-23	Supervisor's initials : CS
-----------------------------------	----------------------------

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.