

HTRF® Research Reagents

HTRF SYK p-Y525/526 kit - 10K pts

Product Number : 64SYKY525PEH Lot Number : 06D

Expiration date : 2025-05-30 Document version : 1

Reagents :	Lot n°
HTRF SYK p-Y525/526 kit - 10K pts d2 antibody	06RA
HTRF SYK p-Y525/526 kit - 10K pts Eu Cryptate antibody	06RA
HTRF SYK p-Y525/526 kit - Ctrl lysate	06RA
HTRF SYK p-Y525/526 kit - modop	1
HTRF P-T prot. - Lysis Buf.3 (4X) 130 mL	14RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	28A
HTRF P-T prot. - Detect. Buf. 50 mL	32RA

Results :

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

Assay characteristics :

	Found value	Specifications Rev 01
Negative	343	/
Control Lysate	3899	/
S/B	11.4	>5

HTRF SYK p-Y525/526 kit - 10K pts - Lot n° : 06D

 Accepted

 Rejected

Quality control date : 2023-01-30	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 USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.