

HTRF® Research Reagents

## HTRF TYK2 p-Y1054/1055 Kit - 10K pts

Product Number : 64TYK2Y10PEH Lot Number : 03A

Expiration date : 15/12/2024 Document version : 1

Reagents :	Lot n°
HTRF TYK2 p-Y1054/1055 Kit - 10K pts d2 antibody	03A
HTRF TYK2 p-Y1054/1055 Kit - 10K pts Eu Cryptate antibody	03A
HTRF TYK2 p-Y1054/1055 Kit - Ctrl Lysate	01RA
HTRF P-T prot. - Lysis Buf.4 (4X) 130 mL	14RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	28B
HTRF P-T prot. - Detect. Buf. 50 mL	32RA

**Results :**

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

**Assay characteristics :**

	Found value	Specifications Rev 01
Negative	389	/
Control Lysate	2138	/
S/B	5.5	>5
	/	/

HTRF TYK2 p-Y1054/1055 Kit - 10K pts - Lot n° : 03A

 Accepted

 Rejected

Quality control date : 2024-04-04	Supervisor's initials : AC
-----------------------------------	----------------------------

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.