

HTRF® Research Reagents

HTRF TYK2 Total Kit - 500 pts

Product Number : 64TYK2TPEG Lot Number : 01E

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

Reagents :	Lot n°
HTRF TYK2 Total Kit - 500 pts d2 antibody	01A
HTRF TYK2 Total Kit - 500 pts Eu Cryptate antibody	01A
HTRF TYK2 Total Kit - Ctrl Lysate	01A
HTRF TYK2 Total Kit - modop	1
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	10E
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	25A
HTRF P-T prot. - Detect. Buf. 2 mL	28C

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	479	/
Control Lysate	3467	/
S/B	7,2	>5
	/	/

HTRF TYK2 Total Kit - 500 pts - Lot n° : 01E

Accepted

Rejected

Quality control date : 2023-04-05	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.