

HTRF® Research Reagents

HTRF STAT3 p-Y705 kit - 10K pts

 Product Number : **62AT3PEH** Lot Number : **22D**

 Expiration date : **14/02/2025** Document version : **1**

Reagents :	Lot n°
HTRF STAT3 p-Y705 kit - 10K pts d2 antibody	22RA
HTRF STAT3 p-Y705 kit - 10K pts Eu Cryptate antibody	22RA
HTRF STAT3 p-Y705 kit - Ctrl lysate	35AR
HTRF P-T prot. - Lysis Buf.1 (4X) 130 mL	34RA
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 : **/**
 Fitting model : **/**
Assay characteristics :

	Found value	Specifications Rev 01
Negative	424	/
Control Lysate	2396	/
S/B	5,7	>5
	/	/

HTRF STAT3 p-Y705 kit - 10K pts - Lot n° : 22D
 Accepted

 Rejected

Quality control date : 2024-01-16	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.