

HTRF® Research Reagents

HTRF STAT3 p-Y705 kit - 96 pts

Product Number : 62AT3PET

Lot Number : 23H

Expiration date : 2025-06-12

Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
HTRF STAT3 p-Y705 kit - Ctrl lysate	35AR
HTRF STAT3 p-Y705 kit - 96 pts d2 antibody	23RA
HTRF STAT3 p-Y705 kit - 96 pts Eu Cryptate antibody	23RA
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	35B
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28B
HTRF P-T prot. - Detect. Buf. 0.5 mL	23RB

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	380	/
Control Lysate	2507	/
S/B	6.6	>5
	/	/

HTRF STAT3 p-Y705 kit - 96 pts - Lot n° : 23H

☒ 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.