

HTRF® Research Reagents

HTRF STAT3 p-Y705 kit - 96 pts

Product Number : 62AT3PET Lot Number : 23J

Expiration date : 2025-06-12 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2413
HTRF STAT3 p-Y705 kit - Ctrl lysate	36A
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	410	/
Control Lysate	3106	/
S/B	7.6	>5
	/	/

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