

HTRF® Research Reagents

HTRF STAT3 total kit - 500 pts

Product Number : 64NT3PEG Lot Number : 05E

Expiration date : 2024-11-20 Document version : 1

Reagents :	Lot n°
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	31F
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	25A
HTRF P-T prot. - Detect. Buf. 2 mL	29A
HTRF STAT3 total kit - 500 pts d2 antibody	05RB
HTRF STAT3 total kit - 500 pts Eu Cryptate antibody	05RB
HTRF STAT3 total kit - Ctrl lysate	08RA
HTRF STAT3 total kit - modop	6

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	409	/
Control Lysate	4117	/
S/B	10,1	>5
/	/	/

HTRF STAT3 total kit - 500 pts - Lot n° : 05E

 Accepted

 Rejected

Quality control date : 2022-08-02	Supervisor's initials : PR
-----------------------------------	----------------------------

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.