

HTRF® Research Reagents

HTRF p70S6K p-T389 kit - 500 pts

Product Number : 64S6KPEG Lot Number : 06F

Expiration date : 2025-12-05 Document version : 1

Reagents :	Lot n°
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	35AR
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28RA
HTRF P-T prot. - Detect. Buf. 2 mL	31RA
HTRF p70S6K p-T389 kit - 500 pts d2 antibody	06RA
HTRF p70S6K p-T389 kit - 500 pts Eu Cryptate antibody	06RA
HTRF p70S6K p-T389 kit - Ctrl lysate	09AR

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	485	/
Control Lysate	3868	/
S/B	8	>5
	/	/

HTRF p70S6K p-T389 kit - 500 pts - Lot n° : 06F

Accepted

Rejected

Quality control date : 2023-10-04	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.