

HTRF® Research Reagents

HTRF SMAD3 p-S423/425 kit - 500 pts

Product Number : 63ADK025PEG Lot Number : 07D

Expiration date : 2026-04-03 Document version : 1

Reagents :	Lot n°
HTRF SMAD3 p-S423/425 kit - Ctrl lysate	13A
HTRF SMAD3 p-S423/425 kit - 500 pts d2 antibody	04F
HTRF SMAD3 p-S423/425 kit - 500 pts Eu Cryptate antibody	07A
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	14B
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28B
HTRF P-T prot. - Detect. Buf. 2 mL	32B

Results :

 Incubation : 4h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : /
 Fitting model : /

Assay characteristics :

	Found value	Specifications Rev 01
Negative	396	/
Control Lysate	4753	/
S/B	12	>5
	/	/

HTRF SMAD3 p-S423/425 kit - 500 pts - Lot n° : 07D

 Accepted

 Rejected

Quality control date : 2024-07-03	Supervisor's initials : PL
-----------------------------------	----------------------------

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.