

HTRF® Research Reagents

HTRF Alpha-SMA kit - 500 pts

 Product Number : **62ASMAPEG** Lot Number : **04T**

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

Reagents :	Lot n°
HTRF Alpha-SMA kit - 500 pts d2 antibody	04D
HTRF Alpha-SMA kit - 500 pts Eu Cryptate antibody	04D
HTRF Alpha-SMA kit - Ctrl lysate	05A
HTRF P-T prot. - Lysis Buf.3 (4X) 2 mL	14A
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28B
HTRF P-T prot. - Detect. Buf. 2 mL	32B

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	432	/
Control Lysate	30365	/
S/B	7.1	>5
	/	/

HTRF Alpha-SMA kit - 500 pts - Lot n° : 04T
 Accepted

 Rejected

Quality control date :	2024-03-14	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.