

## HTRF® Research Reagents

## HTRF SMAD2 p-S465/467 kit - 500 pts

Product Number : 64SMAD2S5PEG Lot Number : 01G

Expiration date : 2027-04-14 Document version : 1

Reagents :	Lot n°
HTRF SMAD2 p-S465/467 kit - 500 pts d2 antibody	01B
HTRF SMAD2 p-S465/467 kit - 500 pts Eu Cryptate antibody	01B
HTRF SMAD2 p-S465/467 kit - Ctrl lysate	04A
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	14B
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	29A
HTRF P-T prot. - Detect. Buf. 2 mL	33A

## 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	500	/
Control Lysate	3688	/
S/B	7,4	>5

HTRF SMAD2 p-S465/467 kit - 500 pts - Lot n° : 01G

☒ Accepted☐ Rejected

Quality control date : 2025-03-04	Supervisor's initials : LM
-----------------------------------	----------------------------

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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.