

HTRF® Research Reagents

HTRF SMAD2 p-S465/467 kit - 500 pts

Product Number : 64SMAD2S5PEG Lot Number : 04A

Expiration date : 2027-11-07 Document version : 1

Reagents :	Lot n°
HTRF SMAD2 p-S465/467 kit - 500 pts d2 antibody	04A
HTRF SMAD2 p-S465/467 kit - 500 pts Eu Cryptate antibody	04A
HTRF SMAD2 p-S465/467 kit - Ctrl lysate	05A
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	15A
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 : EnVision
Fitting model : /

Assay characteristics :

	Found value	Specifications Rev 01
Negative	440	/
Control Lysate	3327	/
S/B	7.6	>5

HTRF SMAD2 p-S465/467 kit - 500 pts - Lot n° : 04A

☒ Accepted

☐ Rejected

Quality control date : 2025-08-05 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 USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.