

HTRF® Research Reagents

## HTRF YAP total kit - 500 pts

Product Number : 64YATPEG Lot Number : 05D

Expiration date : 07/07/2024 Document version : 1

Reagents :	Lot n°
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	12RJ
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	26RA
HTRF P-T prot. - Detect. Buf. 2 mL	29RA
HTRF YAP total kit - 500 pts d2 antibody	05RA
HTRF YAP total kit - 500 pts Eu Cryptate antibody	05RA
HTRF YAP total kit - Ctrl lysate	10RA

**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	400	/
Control Lysate	3547	/
S/B	8,9	>5
	/	/

**HTRF YAP total kit - 500 pts - Lot n° : 05D**
 Accepted

 Rejected

Quality control date : 2021-07-08 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.