

## HTRF® Research Reagents

**HTRF Ubiquitin p-S65 kit - 10K pts**

Product Number : 64UBIS65PEH Lot Number : 04C

Expiration date : 2025-04-05 Document version : 1

Reagents :	Lot n°
HTRF Activation Buffer - 20 mL	04RA
HTRF Ubiquitin p-S65 kit - 10K pts d2 antibody	04RA
HTRF Ubiquitin p-S65 kit - 10K pts Eu Cryptate antibody	04RA
HTRF Ubiquitin p-S65 kit - Ctrl lysate	02RA
HTRF P-T prot. - Lysis Buf.1 (4X) 130 mL	34RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	23D
HTRF P-T prot. - Detect. Buf. 50 mL	32RA

**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	438	/
Control Lysate	7004	/
S/B	16	>5
	/	/

**HTRF Ubiquitin p-S65 kit - 10K pts - Lot n° : 04C**
 Accepted

 Rejected

Quality control date : 2022-11-30 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.