

HTRF® Research Reagents

## HTRF (h) CHK2 p-T68 Kit - 500 pts

Product Number : 64CHK2T8PEG Lot Number : 02B

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

Reagents :	Lot n°
HTRF (h) CHK2 p-T68 Kit - 500 pts d2 antibody	02AR
HTRF (h) CHK2 p-T68 Kit - 500 pts Eu Cryptate antibody	02AR
HTRF (h) CHK2 p-T68 Kit - Ctrl Lysate	02AR
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	35B
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 : Envision  
 Fitting model : /

**Assay characteristics :**

	Found value	Specifications Rev 01
Negative	445	/
Control Lysate	5427	/
S/B	12.2	>5
	/	/

HTRF (h) CHK2 p-T68 Kit - 500 pts - Lot n° : 02B

 Accepted

 Rejected

Quality control date : 2023-11-28 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.