

HTRF® Research Reagents

HTRF PARP cleaved-Asp214 kit - 500 pts

 Product Number : **64PARPEG** Lot Number : **02E**

 Expiration date : **20/03/2025** Document version : **1**

Reagents :	Lot n°
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	14A
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28RA
HTRF P-T prot. - Detect. Buf. 2 mL	32A
HTRF PARP cleaved-Asp214 kit - 500 pts d2 antibody	02RA
HTRF PARP cleaved-Asp214 kit - 500 pts Eu Cryptate antibody	02RA
HTRF PARP cleaved-Asp214 kit - Ctrl lys.	02RA

Results :

 Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : /
 Fitting model : /

Assay characteristics :

	Found value	Specifications Rev 01
Control Lysate	2894	2091-3510 (RATIO)
S/B	6,6	>5

HTRF PARP cleaved-Asp214 kit - 500 pts - Lot n° : 02E

 Accepted

 Rejected

Quality control date : 2024-01-17	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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.