

## HTRF® Research Reagents

**HTRF (h/m) BRD9 Total Kit - 500 pts**

 Product Number : **64BRD9TPEG** Lot Number : **01C**

 Expiration date : **24/02/2025** Document version : **1**

Reagents :	Lot n°
HTRF (h/m) BRD9 Total Kit - 500 pts d2 antibody	01RA
HTRF (h/m) BRD9 Total Kit - 500 pts Eu Cryptate antibody	01RA
HTRF (h/m) BRD9 Total Kit - Ctrl Lysate	01RA
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	14A
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	26RA
HTRF P-T prot. - Detect. Buf. 2 mL	32A

**Results :**

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

	Found value	Specifications Rev 01
<b>Negative</b>	419	/
<b>Control Lysate</b>	5046	/
<b>S/B</b>	12	>5
	/	/

**HTRF (h/m) BRD9 Total Kit - 500 pts - Lot n° : 01C**
 Accepted

 Rejected

Quality control date : <b>2023-03-24</b>	Supervisor's initials : <b>AC</b>
--	-----------------------------------

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.