

## HTRF® Research Reagents

**HTRF (h/m) RAB10 Total Kit - 500 pts**

Product Number : 64RAB10TPEG Lot Number : 01G

Expiration date : 2025-08-01 Document version : 1

Reagents :	Lot n°
HTRF Phospho & Total Activation buffer A - 1.8 mL	05RA
HTRF Phospho & Total Activation buffer B - 1.8 mL	05C
HTRF (h/m) RAB10 Total Kit - 500 pts d2 antibody	01B
HTRF (h/m) RAB10 Total Kit - 500 pts Eu Cryptate antibody	01B
HTRF (h/m) RAB10 Total Kit - Ctrl Lysate	01RA
HTRF P-T prot. - Lysis Buf.3 (4X) 2 mL	14B
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 : Pherastar Plus  
 Fitting model : /

**Assay characteristics :**

	Found value	Specifications Rev 01
Negative	451	/
Control Lysate	3718	/
S/B	8.2	<5
	/	/

HTRF (h/m) RAB10 Total Kit - 500 pts - Lot n° : 01G

 Accepted

 Rejected

Quality control date : 2024-11-26	Supervisor's initials : PR
-----------------------------------	----------------------------

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.