

HTRF® Research Reagents

HTRF cRaf Total Kit - 10K pts

Product Number : 64CRAFTPEH Lot Number : 17A

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

Reagents :	Lot n°
HTRF cRaf Total Kit - 10K pts d2 antibody	17A
HTRF cRaf Total Kit - 10K pts Eu Cryptate antibody	17A
HTRF cRaf Total Kit - Ctrl lysate	15A
HTRF P-T prot. - Lysis Buf.2 (4X) 130 mL	13AR
HTRF P-T prot. - Block. reag.(100X) 2 mL	28B
HTRF P-T prot. - Detect. Buf. 50 mL	32RA

Results :

 Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Envision
 Fitting model : 8

Assay characteristics :

	Found value	Specifications Rev 01
Negative	466	/
Control Lysate	5213	/
S/B	11.2	>5

HTRF cRaf Total Kit - 10K pts - Lot n° : 17A

 Accepted

 Rejected

Quality control date : 2024-07-12	Supervisor's initials : ML
-----------------------------------	----------------------------

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.