

HTRF® Research Reagents

HTRF (h) KRAS Total Kit - 10K pts

Product Number : 64KRASTPEH Lot Number : 05A

Expiration date : 2027-11-07 Document version : 1

Reagents :	Lot n°
HTRF (h) KRAS Total Kit - 10K pts d2 antibody	05A
HTRF (h) KRAS Total Kit - 10K pts Eu Cryptate antibody	05A
HTRF (h) KRAS Total Kit - Ctrl Lysate	03A
HTRF P-T prot. - Lysis Buf.6 (4X) 130 mL	03A
HTRF P-T prot. - Block. reag.(100X) 2 mL	29A
HTRF P-T prot. - Detect. Buf. 50 mL	33A

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	463	/
Control Lysate	4308	/
S/B	9,3	>5

HTRF (h) KRAS Total Kit - 10K pts - Lot n° : 05A

☒ Accepted

☐ Rejected

Quality control date : 2025-05-12	Supervisor's initials : LM
-----------------------------------	----------------------------

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.