

HTRF® Research Reagents

HTRF IRAK4 total kit - 500 pts

Product Number : 63ADK108PEG Lot Number : 09D

Expiration date : 2024-09-14 Document version : 1

Reagents :	Lot n°
HTRF IRAK4 total kit - 500 pts d2 antibody	09A
HTRF IRAK4 total kit - 500 pts Eu Cryptate antibody	09A
HTRF IRAK4 total kit - Ctrl lysate	09RA
HTRF IRAK4 total kit - modop	5
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	31F
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	26A
HTRF P-T prot. - Detect. Buf. 2 mL	28C

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	433	/
Control Lysate	4202	/
S/B	9,7	>5
/	/	/

HTRF IRAK4 total kit - 500 pts - Lot n° : 09D

 Accepted

 Rejected

Quality control date : 2022-06-29	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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.