

HTRF® Research Reagents

HTRF Human ApoB Kit - 500 pts

Product Number : 64APOBPEG Lot Number : 01A

Expiration date : 2025-06-16 Document version : 1

Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20RA
HTRF Detect. Buf. 3 - 7 mL	30AR
HTRF Detect. Buf. 3 - 7 mL	30AR
HTRF Human ApoB Kit - Std	01A
HTRF Human ApoB Kit - 500 pts d2 antibody	01A
HTRF Human ApoB Kit - 500 pts Eu Cryptate antibody	01A

Results :

Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : /
 Fitting model : nonlinear regression, sigmoïdal dose response (variable slope)

Assay characteristics :

	mean - 2SD	mean + 2SD	Soit +/-
delta F %			
10.3	7	16	20%
27.90	14	41	25%
75.3	55	85	20%
203	43	325	38%
549	278	454	20%
1481	1658	2407	20%
4000	3998	4727	20%
LOQ	11	51	32%

HTRF Human ApoB Kit - 500 pts - Lot n° : 01A

Accepted

Rejected

Quality control date : 2024-03-06 Supervisor's initials : AC

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.