

HTRF® Research Reagents

HTRF Aldosterone kit - 500 pts

Product Number : 64ALDPEG Lot Number : 09C

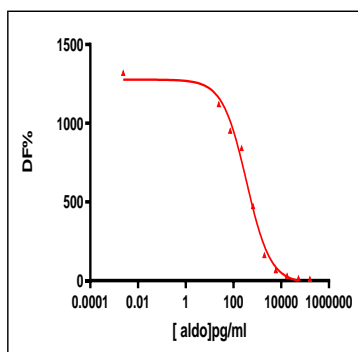
Expiration date : 16/06/2025 Document version : 1

Reagents :	Lot n°
Aldosterone kit standard	05A
Aldosterone kit -500 tests Eu Cryptate antibody	07C
Aldosterone kit - 500 tests XL665 reagent	09A
Aldosterone kit - modop	4
Diluent 1 - 20 mL	19A
Detection buffer 3 - 7 mL	20A

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

ALDO	Conc. (ng/ml)	DF %
Std 0	0	1319%
Std 1	0,024	1121%
Std 2	0,073	952%
Std 3	0,22	842%
Std 4	0,65	473%
Std 5	1,9	162%
Std 6	5,9	66%
Std 7	17,7	30%
Std 8	53,3	15%
Std 9	160	10%



	Found value	Specifications Rev 01
DF max %	1319	1000 – 1700 %
EC 50	350 pg/ml	250 – 350 pg/ml
detection limite	6 pg/ml	≤ 25 pg/ml

HTRF Aldostérone kit - 500 pts - Lot n° : 09C

Accepted

Rejected

Quality control date :	2023-09-15	Supervisor's initials :	FMV
------------------------	------------	-------------------------	-----

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