

HTRF® Research Reagents

## HTRF Aldosterone kit - 500 pts

Product Number : 64ALDPEG Lot Number : 10C

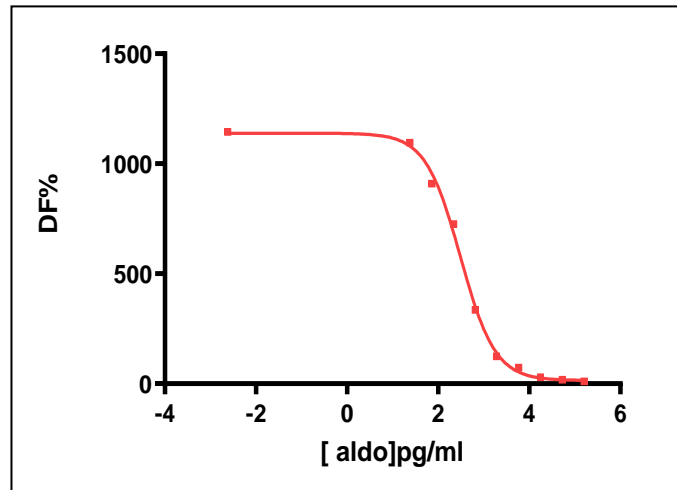
Expiration date : 2025-06-16 Document version : 1

Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20RA
HTRF Detect. Buf. 3 - 7 mL	30AR
HTRF Detect. Buf. 3 - 7 mL	30AR
HTRF Aldosterone kit - Std	07A
HTRF Aldosterone kit - 500 pts Eu Cryptate antibody	07D
HTRF Aldosterone kit - 500 pts XL665 reagent	10B

**Results :**

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

ALDO	Conc. (ng/ml)	DF %
Std 0	0	1144%
Std 1	0,024	1096%
Std 2	0,073	909%
Std 3	0,22	726%
Std 4	0,65	336%
Std 5	1,9	124%
Std 6	5,9	73%
Std 7	17,7	29%
Std 8	53,3	18%
Std 9	160	10%



	Found value	Specifications Rev 01
DF max %	1144	1000 – 1700 %
EC 50	309 pg/ml	250 – 350 pg/ml
detection limite	23	≤ 25 pg/ml

HTRF Aldosterone kit - 500 pts - Lot n° : 10C

Accepted

Rejected

Quality control date : 2024-07-18	Supervisor's initials : PL
-----------------------------------	----------------------------

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.