

Certificate of Analysis

PROGEST-RIA PROGESTERONE RIA

Lot Nr : 213306

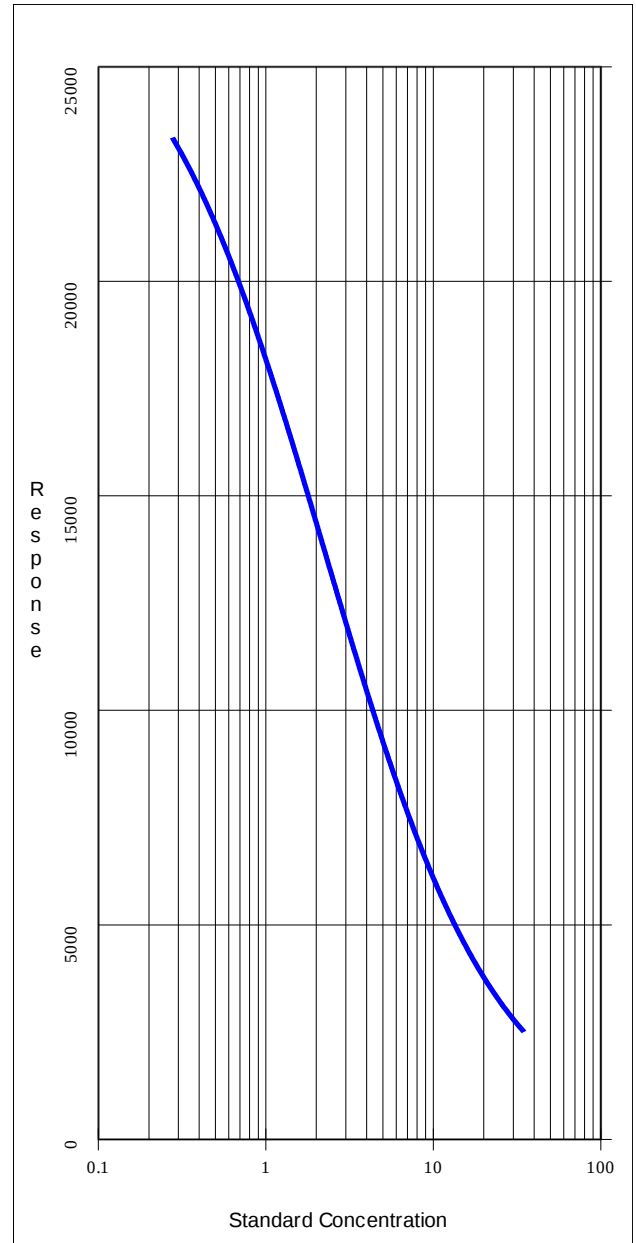
Expiry Date : 29/10/2021

Tested composition :

<i>Description</i>	<i>Lot nr.</i>	<i>Expiry Date</i>
TRAC	21H09/02	29/10/2021
TUBES	2108-3017/02	01/08/2024
STD 0	1902-2404/04	19/04/2024
C1	1812-1695/06	17/08/2023
C2	1812-1697/03	17/08/2023
STD	1906-3989/02	17/10/2024
WASH-SOL	2101-1059/04	26/12/2025

Results:

<i>Parameter</i>	<i>Response</i>	<i>B/T</i>	<i>B/B0</i>	<i>Conc</i>	<i>Range</i>
	<i>Cpm</i>	<i>%</i>	<i>%</i>	<i>ng/ml</i>	<i>ng/ml</i>
TOT	62976.0				
STD0 1902-2404/04	24706.5	39.23			
STD1 1906-3989/02	23642.0	37.54	95.69	0.25	
STD2 1906-3989/02	18854.0	29.94	76.31	0.85	
STD3 1906-3989/02	11795.0	18.73	47.74	3.32	
STD4 1906-3989/02	6496.7	10.32	26.30	8.74	
STD5 1906-3989/02	4383.0	6.96	17.74	16.3	
STD6 1906-3989/02	2513.3	3.99	10.17	34.8	
C1 1812-1695 06	18263.7	29.00	73.92	0.98	1.12 ± 0.48
C2 1812-1697 03	6505.7	10.33	26.33	9.03	8.43 ± 2.52



Only the figures before the slash should be taken into account for the batch identification.

Verified by : **Nicolas Grégoire**

Edition date : **10/08/2021**

ACCEPTED

This document has been produced electronically and is valid without a signature.