

revvity

Biomedical Waste Disposal Record

RHS\_F\_GN\_028


v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

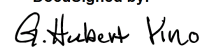
Month/ Year: April 2026

S. No.	Date of Generation	Quantity of BMW Generated (in KG) Colour Coding and Category					Date of collection by Agency	Time (in AM/ PM)	Name of Waste Collector	Name of Staff
		Yellow	Red	White	Blue	Total				
1.	31-03-2026	-	42	-	4	46	01-04-2026	10:55am	Saravanan	Jintu Das
2.	02-04-2026	-	16	-	-	16	03-04-2026	10:30am	Saravanan	Rajkiran
3.	04-04-2026	-	13	-	-	13	06-04-2026	10:40am	Saravanan	Rajkiran
4.	07-04-2026	-	38	-	-	38	08-04-2026	10:30am	Saravanan	Rajkiran
5.	09-04-2026	-	14	-	-	14	10-04-2026	10.30am	Saravanan	Rajkiran
6.	11-04-2026	-	11	-	2	13	13-04-2026	11:00am	Saravanan	Rajkiran
7.	13-04-2026	-	16	-	-	16	15-04-2026	10:30am	Saravanan	Rajkiran
8.	16-04-2026	-	17	-	3	20	17-04-2026	10:30am	Saravanan	Rajkiran
9.	18-04-2026	-	15	-	-	15	20-04-2026	10:10am	Madhan	Rajkiran
10.	21-04-2026	-	32	-	-	32	22-04-2026	10:30am	Saravanan	Rajkiran
11.	22-04-2026	-	17	-	-	17	24-04-2026	10:40am	Saravanan	Rajkiran
12.	25-04-2026	-	18	-	5	23	27-04-2026	10:10am	Saravanan	Rajkiran
13.	28-04-2026	-	54	-	-	54	29-04-2026	10:05am	Saravanan	Rajkiran
14.										
	<b>TOTAL (Kgs)</b>	-	<b>303</b>	-	<b>14</b>	<b>317</b>				

DocuSigned by:  
  
 DC2050D1B1DE4DF... (Sign & Date)

Checked by

04 May 2026 | 3:33 AM EDT

DocuSigned by:  
  
 02CF50F7F1A44D8... (Sign & Date)

Verified by

04 May 2026 | 4:15 AM EDT