

revvity

Biomedical Waste Disposal Record

RHS\_F\_GN\_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: May 2024

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.	02-05-2024	-	20	-	-	20	03-05-2024	10:00am	Raghu	Deepak
2.	04-05-2024	7	16	-	-	23	06-05-2024	10:00am	Sreeajith	Suman Banik
3.	07-05-2024	-	35	-	10	45	08-05-2024	10:30am	Saravanan	Suman Banik
4.	09-05-2024	-	15	-	2	17	10-05-2024	09:20am	Raghu	Suman Banik
5.	11-05-2024	9	24	-	-	33	13-05-2024	10.05am	Muthu Kumar	Deepak
6.	14-05-2024	9	16	-	-	25	15-05-2024	10:20am	Muthu Kumar	Deepak
7.	16-05-2024	7	14	-	3	24	17-05-2024	09:50am	Muthu Kumar	Deepak
8.	18-05-2024	-	18	-	-	18	20-05-2024	09:59am	Muthu Kumar	Suman Banik
9.	21-05-2024	6	13	-	-	19	22-05-2024	11:00am	Muthu Kumar	Suman Banik
10.	23-05-2024	7	14	-	-	21	24-05-2024	11:05am	Muthu Kumar	Suman Banik
11.	25-05-2024	12	31	-	-	43	27-05-2024	10:15am	Saravanan	Deepak
12.	28-05-2024	11	47	-	6	64	29-05-2024	10:00am	Saravanan	Deepak
13.	30-05-2024	6	28	-	-	34	31-05-2024	09:50am	Saravanan	Deepak
	<b>TOTAL (Kgs)</b>	<b>74</b>	<b>291</b>		<b>21</b>	<b>386</b>				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF... (Sign &amp; Date)

Checked by

03 June 2024 | 8:53 AM EDT

DocuSigned by:

G. Hubert Vino

02CE50E7E1A44D8 (Sign &amp; Date)

Verified by

03 June 2024 | 1:31 PM EDT