

revvity

Biomedical Waste Disposal Record

RHS\_F\_GN\_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: September 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.	31-08-2024	-	14	-	5	19	02-09-2024	10:05am	Gopinath	Deepak
2.	03-09-2024	2	41	-	5	48	04-09-2024	10:20am	Saravanan	Deepak
3.	05-09-2024	1	17	-	3	21	06-09-2024	11:05am	Saravanan	Deepak
4.	07-09-2024	1	17	-	10	28	09-09-2024	10:30am	Saravanan	Suman Banik
5.	10-09-2024	-	19	-	-	19	11-09-2024	10.25am	Saravanan	Suman Banik
6.	12-09-2024	1	24	-	2	27	13-09-2024	10:00am	Saravanan	Suman Banik
7.	14-09-2024	-	42	-	-	42	16-09-2024	10:18am	Saravanan	Deepak
8.	17-09-2024	-	21	-	4	25	18-09-2024	10:30am	Sivaraj	Deepak
9.	19-09-2024	-	31	-	-	31	20-09-2024	10:10am	Saravanan	Deepak
10.	21-09-2024	1	27	-	-	28	23-09-2024	10:07am	Saravanan	Suman Banik
11.	24-09-2024	-	17	-	2	19	25-09-2024	10:10am	Saravanan	Deepak
12.	27-09-2024	-	16	-	-	16	27-09-2024	10:40am	Saravanan	Deepak
13.	28-09-2024	-	34	-	-	34	30-09-2024	10:00am	Abin	Deepak
14.										
	<b>TOTAL (Kgs)</b>	<b>6</b>	<b>320</b>		<b>31</b>	<b>357</b>				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF...

(Sign &amp; Date)

Checked by

30 September 2024 | 3:55 AM EDT

DocuSigned by:

G. Hubert Kuo

02CF50F7F1A44D8...

(Sign &amp; Date)

Verified by

30 September 2024 | 9:28 AM EDT