


revvity	<i>Biomedical Waste Disposal Record</i>	RHS_F_GN_028	v.1
Effective Date: 01/January/2024		Approval Date: 18/February/2019	

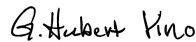
Month/ Year: June 2024

S. No.	Date of Generation	Quantity of BMW Generated (in KG)					Date of collection by Agency	Time (in AM/ PM)	Name of Waste Collector	Name of Staff
		Colour Coding and Category								
		Yellow	Red	White	Blue	Total				
1.	01-06-2024	4	12	-	-	16	03-06-2024	10:17am	Saravanan	Suman Banik
2.	04-06-2024	5	11	-	-	16	05-06-2024	10:56am	Saravanan	Suman Banik
3.	06-06-2024	3	11	-	-	14	07-06-2024	10:41am	Saravanan	Suman Banik
4.	08-06-2024	4	11	-	-	15	10-06-2024	12:05pm	Saravanan	Deepak
5.	11-06-2024	6	18	-	-	24	12-06-2024	10.45am	Saravanan	Deepak
6.	13-06-2024	12	31	-	4	47	14-06-2024	10:00am	Saravanan	Deepak
7.	15-06-2024	6	17	-	-	23	17-06-2024	10:00am	Saravanan	Suman Banik
8.	18-06-2024	6	28	-	-	34	19-06-2024	10:15am	Saravanan	Suman Banik
9.	20-06-2024	5	17	-	-	22	21-06-2024	09:55am	Saravanan	Suman Banik
10.	22-06-2024	14	37	-	-	51	24-06-2024	10:20am	Saravanan	Deepak
11.	25-06-2024	5	17	-	-	22	26-06-2024	10:15am	Saravanan	Deepak
12.	27-06-2024	13	21	-	-	34	28-06-2024	10:30am	Saravanan	Deepak
	TOTAL (Kgs)	83	231		4	318				

DocuSigned by:

 DC2050D1B1DE4DF... (Sign & Date)

Checked by

01 July 2024 | 4:25 AM EDT

DocuSigned by:

 02CF50F7F1A44D8... (Sign & Date)

Verified by

01 July 2024 | 1:56 PM EDT