

revvity

Biomedical Waste Disposal Record

RHS_F_GN_028


v.1

Effective Date: 01/January/2024

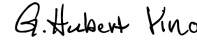
Approval Date: 18/February/2019

Month/ Year: July 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.	29-06-2024	4	15	-	-	19	01-07-2024	10:04am	Arun	Suman Banik
2.	02-07-2024	13	31	-	-	44	03-07-2024	10:38am	Saravanan	Suman Banik
3.	04-07-2024	13	31	-	-	44	05-07-2024	10:55am	Saravanan	Suman Banik
4.	06-07-2024	4	17	-	-	21	08-07-2024	10:35pm	Saravanan	Deepak
5.	09-07-2024	6	17	-	4	27	10-07-2024	10.20am	Saravanan	Deepak
6.	11-07-2024	6	13	-	-	19	12-07-2024	10:18am	Saravanan	Deepak
7.	13-07-2024	7	14	-	3	24	15-07-2024	10:10am	Saravanan	Suman Banik
8.	16-07-2024	-	15	-	-	15	17-07-2024	10:00am	Saravanan	Suman Banik
9.	18-07-2024	11	31	-	-	42	19-07-2024	11:19am	Saravanan	Suman Banik
10.	20-07-2024	12	31	-	-	43	22-07-2024	10:15am	Saravanan	Deepak
11.	23-07-2024	5	13	-	-	18	24-07-2024	10:00am	Saravanan	Deepak
12.	25-07-2024	-	18	-	4	22	26-07-2024	10:20am	Saravanan	Deepak
13.	27-07-2024	5	15	-	-	20	29-07-2024	10:14am	Abin	Suman Banik
14.	30-07-2024	1	12	-	-	13	31-07-2024	10:50am	Saravanan	Suman Banik
	TOTAL (Kgs)	87	273		11	371				

DocuSigned by:

 DC2050D1B1DE4DF... (Sign & Date)

Checked by
 31 July 2024 | 4:35 AM EDT

DocuSigned by:

 02CF50F7F1A44D8... (Sign & Date)

Verified by
 31 July 2024 | 5:13 AM EDT