
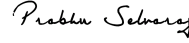


<b>revvity</b>	<i>Biomedical Waste Disposal Record</i>	<b>RHS_F_GN_028</b>	<b>v.1</b>
Effective Date: 01/January/2024		Approval Date: 18/February/2019	

**Month/ Year: April 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.	30-03-2024	8	43	-	-	51	01-04-2024	09:40 am	Saravanan	Deepak
2.	02-04-2024	6	16	-	4	26	03-04-2024	10:15 am	Abin	Deepak
3.	04-04-2024	6	10	-	2	18	05-04-2024	09:30 am	Abin	Deepak
4.	06-04-2024	3	12	-	2	17	08-04-2024	09:40 am	Abin	Suman Banik
5.	09-04-2024	8	27	-	4	39	10-04-2024	09:59 am	Saravanan	Suman Banik
6.	11-04-2024	8	13	-	-	21	12-04-2024	10:45 am	Saravanan	Suman Banik
7.	13-04-2024	6	18	-	-	24	15-04-2024	09:40 am	Saravanan	Deepak
8.	16-04-2024	5	28	-	-	33	17-04-2024	10:05 am	Saravanan	Deepak
9.	18-04-2024	7	29	-	-	36	22-04-2024	09:45 am	Saravanan	Suman Banik
10.	23-04-2024	3	20	-	-	23	24-04-2024	09:35 am	Abin	Suman Banik
11.	25-04-2024	7	32	-	-	39	26-04-2024	09:50 am	Saravanan	Suman Banik
12.	27-04-2024	6	21	-	-	27	29-04-2024	09:58 am	Saravanan	Deepak
	<b>TOTAL (Kgs)</b>	<b>73</b>	<b>269</b>	<b>-</b>	<b>12</b>	<b>354</b>				

DocuSigned by:  
  
 CSBDD64A5DF1495...  
 08 May 2024 | 5:02 AM EDT  
 (Sign & Date)  
**Checked by QA**

DocuSigned by:  
  
 3849F7FB50A0499...  
 08 May 2024 | 6:37 AM EDT  
 (Sign & Date)  
**Verified by QM**