

<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: August 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.	01-08-2024	3	26	-	-	29	02-08-2024	10:30am	Saravanan	Suman Banik
2.	03-08-2024	-	20	-	5	25	05-08-2024	10:35am	Raghu	Deepak
3.	06-08-2024	1	15	-	-	16	07-08-2024	10:00am	Saravanan	Deepak
4.	08-08-2024	-	29	-	-	29	09-08-2024	10:54pm	Saravanan	Deepak
5.	10-08-2024	-	15	-	-	15	12-08-2024	10.30am	Saravanan	Suman Banik
6.	13-08-2024	-	17	-	-	17	14-08-2024	10:40am	Saravanan	Suman Banik
7.	15-08-2024	-	16	-	-	16	16-08-2024	10:25am	Saravanan	Suman Banik
8.	17-08-2024	1	52	-	-	53	19-08-2024	10:10am	Saravanan	Deepak
9.	20-08-2024	-	14	-	-	14	21-08-2024	10:20am	Saravanan	Deepak
10.	22-08-2024	-	46	-	5	51	23-08-2024	10:50am	Saravanan	Deepak
11.	24-08-2024	-	18	-	-	18	26-08-2024	09:29am	Abin	Suman Banik
12.	27-08-2024	-	8	-	-	8	28-08-2024	09:57am	Raghu	Suman Banik
13.	29-08-2024	-	63	-	-	63	30-08-2024	14:06am	Gopalan	Suman Banik
14.										
	<b>TOTAL (Kgs)</b>	<b>5</b>	<b>339</b>		<b>10</b>	<b>354</b>				

DocuSigned by:  
*Baranidaran S*  
DC2050D1B1DE4DF... (Sign & Date)

**Checked by**

31 August 2024 | 6:29 AM EDT

DocuSigned by:  
*G. Hubert Kuro*  
02CF50F7F1A44D8... (Sign & Date)

**Verified by**

31 August 2024 | 6:43 AM EDT