


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

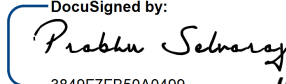
Month/ Year: March 2024

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.	01-03-2024		20			20	01-03-2024	12:40pm	Raghu	Deepak
2.	02-03-2024	7	46		4	57	04-03-2024	10:00am	Saravanan	Deepak
3.	05-03-2024	11	21			32	06-03-2024	10:45am	Saravanan	Deepak
4.	07-03-2024	3	15		4	22	08-03-2024	10:00am	Saravanan	Deepak
5.	09-03-2024	2	14			16	11-03-2024	10.08am	Saravanan	Suman Banik
6.	12-03-2024	7	41			48	13-03-2024	10:10am	Saravanan	Suman Banik
7.	14-03-2024	6	28			34	15-03-2024	09:45am	Saravanan	Suman Banik
8.	16-03-2024	3	17		2	22	18-03-2024	10:30am	Saravanan	Deepak
9.	19-03-2024	8	48		4	60	20-03-2024	10:05am	Saravanan	Deepak
10.	21-03-2024	3	17		4	24	22-03-2024	10:00am	Saravanan	Deepak
11.	23-03-2024	4	14			18	25-03-2024	09:40am	Saravanan	Suman Banik
12.	26-03-2024	5	29			34	27-03-2024	10:10am	Saravanan	Suman Banik
13.	28-03-2024	9	18			27	29-03-2024	09:28am	Abin	Suman Banik
	TOTAL (Kgs)	68	328		18	414				

DocuSigned by:

 A09C56B9BCB349E... (Sign & Date)

Checked by QA

01 April 2024 | 5:27 AM EDT

DocuSigned by:

 3849F7FB50A0499... (Sign & Date)

Verified by QM

04 April 2024 | 1:47 AM EDT