
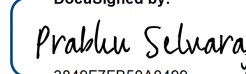


revvity	<i>Biomedical Waste Disposal Record</i>	RHS_F_GN_028	v.1
Effective Date: 01/January/2024	Approval Date: 18/February/2019		

Month/ Year: February 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.	01-02-2024	6	24			30	02-02-2024	10:24am	Saravanan	Suman Banik
2.	03-02-2024	5	14			19	05-02-2024	09:45am	Saravanan	Suman Banik
3.	06-02-2024	6	37			43	07-02-2024	10:26am	Saravanan	Suman Banik
4.	08-02-2024	6	21		4	31	09-02-2024	10:30am	Saravanan	Suman Banik
5.	10-02-2024		20			20	11-02-2024	11.40am	Raghu	Deepak
6.	13-02-2024	8	18			26	14-02-2024	09:40am	Arun	Deepak
7.	15-02-2024	16	47			63	16-02-2024	10:45am	Saravanan	Deepak
8.	17-02-2024		20			20	19-02-2024	12:25am	Raghu	Suman Banik
9.	20-02-2024	8	17		4	29	21-02-2024	10:35am	Saravanan	Suman Banik
10.	22-02-2024	11	27			38	23-02-2024	10:13am	Saravanan	Suman Banik
11.	24-02-2024	11	32			43	26-02-2024	10:00am	Saravanan	Deepak
12.	27-02-2024	7	31			38	28-02-2024	10:20am	Saravanan	Deepak
13.										
	TOTAL (Kgs)	84	308		8	400				

DocuSigned by:

 A09C56B9BCB349E... (Sign & Date)
Checked by QA
 07 March 2024 | 6:42 AM EST

DocuSigned by:

 3849F7FB50A0499... (Sign & Date)
Verified by QM
 20 March 2024 | 3:13 AM EDT