

revvity

Biomedical Waste Disposal Record

RHS_F_GN_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: December 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.	30-11-2024	-	18	-	-	18	02-12-2024	10:15am	Saravanan	Deepak
2.	03-12-2024	-	48	-	3	51	04-12-2024	11:39am	Saravanan	Deepak
3.	05-12-2024	-	16	-	-	16	06-12-2024	10:18am	Saravanan	Deepak
4.	07-12-2024	-	29	-	-	29	09-12-2024	10:16am	Saravanan	Deepak
5.	09-12-2024	-	17	-	-	17	11-12-2024	10.51am	Saravanan	Deepak
6.	12-12-2024	-	16	-	3	19	13-12-2024	11:16am	Saravanan	Deepak
7.	14-12-2024	-	17	-	-	17	16-12-2024	10:46am	Saravanan	Deepak
8.	17-12-2024	-	24	-	-	24	18-12-2024	10:30am	Saravanan	Deepak
9.	19-12-2024	-	27	-	-	27	20-12-2024	10:40am	Saravanan	Deepak
10.	21-12-2024	-	16	-	2	18	23-12-2024	10:15am	Saravanan	Deepak
11.	26-12-2024	-	15	-	-	15	27-12-2024	09:50am	Ragu	Deepak
12.	28-12-2024	-	28	-	-	28	30-12-2024	10:05am	Saravanan	Deepak
13.										
14.										
	TOTAL (Kgs)	-	271	-	8	279				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF...

(Sign & Date)

Checked by

30 December 2024 | 3:01 AM EST

DocuSigned by:

G. Hubert Vino

02CF50F7F1A44D8...

(Sign & Date)

Verified by

30 December 2024 | 5:17 AM EST