

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: January 2025

S. No.	Date of Generation	Quantity of BMW Generated (in KG)					Date of collection by Agency	Time (in AM/ PM)	Name of Waste Collector	Name of Staff
		Colour Coding and Category								
		Yellow	Red	White	Blue	Total				
1.	02-01-2025	-	28	-	-	28	03-01-2025	10:00am	Saravanan	Deepak
2.	04-01-2025	-	27	-	-	27	06-01-2025	10:50am	Saravanan	Deepak
3.	07-01-2025	3	15	-	5	23	08-01-2025	10:30am	Abin	Deepak
4.	09-01-2025	-	32	-	-	32	10-01-2025	10:20am	Abin	Deepak
5.	11-01-2025	-	21	-	3	24	13-01-2025	10.25am	Saravanan	Deepak
6.	16-01-2025	-	23	-	-	23	17-01-2025	10:20am	Saravanan	Deepak
7.	18-01-2025	-	24	-	-	24	20-01-2025	10:22am	Saravanan	Deepak
8.	21-01-2025	-	31	-	-	31	22-01-2025	10:40am	Saravanan	Deepak
9.	23-01-2025	-	17	-	-	17	24-01-2025	10:20am	Saravanan	Deepak
10.	25-01-2025	-	15	-	-	15	27-01-2025	10:20am	Saravanan	Deepak
11.	28-01-2025	-	32	-	-	32	29-01-2025	10:25am	Saravanan	Deepak
12.	31-01-2025	1	41	-	5	27	31-01-2025	10:20am	Saravanan	Deepak
13.										
14.										
	TOTAL (Kgs)	4	306	-	13	323				

DocuSigned by:
Baranidaran S
DC2050D1B1DE4DF...

(Sign & Date)

Checked by

01 February 2025 | 12:18 AM EST

DocuSigned by:
G. Hubert Kuo
02CF50F7F1A44D8...

(Sign & Date)

Verified by

01 February 2025 | 12:46 AM EST