

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 2025

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.	01-02-2025	-	16	-	-	16	03-02-2025	10:45am	Saravanan	Deepak
2.	04-02-2025	-	16	-	-	16	05-02-2025	10:40am	Saravanan	Deepak
3.	06-02-2025	-	24	-	3	27	07-02-2025	01:55pm	Saravanan	Deepak
4.	08-02-2025	-	16	-	-	16	10-02-2025	10:30am	Saravanan	Moorthy
5.	11-02-2025	-	16	-	2	18	12-02-2025	10.15am	Saravanan	Deepak
6.	13-02-2025	-	31	-	1	32	14-02-2025	10:40am	Saravanan	Deepak
7.	15-02-2025	-	36	-	-	36	17-02-2025	10:56am	Saravanan	Deepak
8.	18-02-2025	-	17	-	-	17	19-02-2025	10:40am	Saravanan	Deepak
9.	20-02-2025	-	17	-	-	17	21-02-2025	10:18am	Saravanan	Deepak
10.	22-02-2025	-	32	-	2	34	24-02-2025	10:10am	Saravanan	Deepak
11.	25-02-2025	-	16	-	-	16	26-02-2025	10:30am	Saravanan	Deepak
12.	27-02-2025	-	15	-	-	15	28-02-2025	10:29am	Saravanan	Deepak
13.										
14.										
	TOTAL (Kgs)	-	252	-	8	260				

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

Checked by 28 February 2025 | 12:15 AM EST

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

Verified by 28 February 2025 | 1:32 AM EST