

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: November 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.	03-11-2024	-	26	-	-	26	04-11-2024	10:30am	Saravanan	Deepak
2.	05-11-2024	-	32	-	4	36	06-11-2024	10:20am	Saravanan	Deepak
3.	07-11-2024	-	18	-	1	19	08-11-2024	11:05am	Saravanan	Deepak
4.	09-11-2024	-	24	-	-	24	11-11-2024	10:30am	Saravanan	Deepak
5.	12-11-2024	-	21	-	-	21	13-11-2024	10.15am	Saravanan	Deepak
6.	14-11-2024	-	32	-	2	34	15-11-2024	10:20am	Saravanan	Deepak
7.	16-11-2024	-	24	-	3	27	18-11-2024	10:00am	Saravanan	Deepak
8.	18-11-2024	-	23	-	-	23	20-11-2024	10:28am	Saravanan	Deepak
9.	21-11-2024	-	19	-	-	19	22-11-2024	10:00am	Saravanan	Deepak
10.	24-11-2024	-	21	-	-	21	25-11-2024	10:39am	Saravanan	Moorthy
11.	26-11-2024	-	14	-	3	17	27-11-2024	10:30am	Saravanan	Deepak
12.	28-11-2024	-	17	-	-	17	29-11-2024	10:15am	Saravanan	Deepak
13.										
14.										
	TOTAL (Kgs)	-	271	-	13	284				

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

Checked by 30 November 2024 | 3:37 AM EST

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

Verified by 30 November 2024 | 4:34 AM EST