

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

Month/ Year: April 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.	01-04-2025	-	16	-	-	16	02-04-2025	10:40am	Saravanan	Deepak
2.	03-04-2025	-	17	-	-	17	04-04-2025	10:20am	Saravanan	Deepak
3.	05-04-2025	-	21	-	-	21	07-04-2025	10:43am	Saravanan	Deepak
4.	08-04-2025	-	23	-	3	27	09-04-2025	11:03am	Saravanan	Deepak
5.	10-04-2025	1	24	-	-	25	11-04-2025	10.10am	Saravanan	Deepak
6.	15-04-2025	-	28	-	2	30	16-04-2025	10:30am	Saravanan	Deepak
7.	17-04-2025	-	43.6	-	-	43.6	18-04-2025	10:00am	Ragu	Deepak
8.	19-04-2025	-	18	-	-	18	21-04-2025	10:24am	Saravanan	Moorthy
9.	22-04-2025	2	54	-	-	56	23-04-2025	10:33am	Saravanan	Moorthy
10.	24-04-2025	-	21	-	-	21	25-04-2025	10:40am	Saravanan	Moorthy
11.	26-04-2025	-	16	-	5	21	28-04-2025	10:10am	Saravanan	Deepak
12.	29-04-2025	-	21	-	-	21	30-04-2025	10:40am	Saravanan	Deepak
13.										
14.										
	TOTAL (Kgs)	3	303.6	-	12	316.6				

DocuSigned by:
Baranidaran S
DC2050D1B1DE4DF...

(Sign & Date)

Checked by
30 April 2025 | 1:47 AM EDT

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

(Sign & Date)

Verified by
30 April 2025 | 9:04 AM EDT