



Biomedical Waste Disposal Record

RHS_F_GN_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: July 2026

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.	30-06-2026	-	56	-	-	56	01-07-2026	11:30am	Saravanan	Rajkiran
2.	02-07-2026	-	26	-	-	26	03-07-2026	10:40am	Saravanan	Rajkiran
3.	04-07-2026	-	39	-	3	42	06-07-2026	11:00am	Saravanan	Jintu
4.	07-07-2026	-	24	-	-	24	08-07-2026	10:50am	Saravanan	Jintu
5.	09-07-2026	-	26	-	-	26	10-07-2026	10.20am	Saravanan	Jintu
6.	11-07-2026	-	21	-	3	24	13-07-2026	10:05am	Saravanan	Rajkiran
7.	14-07-2026	-	31	-	-	31	15-07-2026	10:30am	Saravanan	Rajkiran
8.	16-07-2026	-	30	-	-	30	17-07-2026	10:00am	Prakash	Rajkiran
9.	18-07-2026	-	42	-	4	46	20-07-2026	11:05am	Saravanan	Jintu
10.	21-07-2026	-	31	-	-	31	22-07-2026	11:25am	Saravanan	Jintu
11.	23-07-2026	-	19	-	-	19	24-07-2026	10:20am	Saravanan	Jintu
12.	25-07-2026	-	28	-	3	31	27-07-2026	11:00am	Saravanan	Rajkiran
13.	28-07-2026	-	31	-	-	31	29-07-2026	10:45am	Saravanan	Rajkiran
14.	30-07-2026	-	17	-	-	17	31-07-2026	11:00am	Saravanan	Rajkiran
	TOTAL (Kgs)	-	421	-	13	434				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF...

(Sign & Date)

Checked by

01 August 2026 | 8:24 AM EDT

DocuSigned by:

G. Hubert Yno

02CF50F7F1A44D8...

(Sign & Date)

Verified by

01 August 2026 | 9:50 AM EDT