



Biomedical Waste Disposal Record

RHS_F_GN_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: June 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.	30-05-2026	-	28	-	-	28	01-06-2026	10:35am	Saravanan	Rajkiran
2.	02-06-2026	-	24	-	2	26	03-06-2026	10:30am	Saravanan	Rajkiran
3.	04-06-2026	-	42	-	-	42	05-06-2026	10:40am	Saravanan	Rajkiran
4.	06-06-2026	-	38	-	4	42	08-06-2026	11:05am	Saravanan	Jintu
5.	09-06-2026	-	48	-	-	48	10-06-2026	11:15am	Saravanan	Jintu
6.	11-06-2026	-	41	-	-	41	12-06-2026	11:15am	Saravanan	Jintu
7.	13-06-2026	-	52	-	-	52	15-06-2026	10:30am	Saravanan	Rajkiran
8.	16-06-2026	-	41	-	-	41	17-06-2026	10:50am	Saravanan	Rajkiran
9.	18-06-2026	-	37	-	-	37	19-06-2026	10:35am	Saravanan	Rajkiran
10.	20-06-2026	-	42	-	5	47	22-06-2026	11:10am	Saravanan	Rajkiran
11.	23-06-2026	-	19	-	-	19	24-06-2026	10:45am	Saravanan	Jintu
12.	25-06-2026	-	41	-	-	41	26-06-2026	10:46am	Saravanan	Jintu
13.	27-06-2026	-	48	-	4	52	29-06-2026	10:45am	Saravanan	Rajkiran
14.										
	TOTAL (Kgs)	1	501	-	15	516				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF...

(Sign & Date)

Checked by

01 July 2026 | 2:13 AM EDT

DocuSigned by:

G. Hubert Vino

02CF50F7F1A44D8...

(Sign & Date)

Verified by

01 July 2026 | 7:35 AM EDT