



Biomedical Waste Disposal Record

RHS_F_GN_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: May 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.	02-05-2026	-	18	-	2	20	04-05-2026	10:30am	Saravanan	Rajkiran
2.	05-05-2026	-	13	-	-	13	06-05-2026	10:15am	Saravanan	Rajkiran
3.	07-05-2026	-	21	-	-	21	08-05-2026	10:40am	Saravanan	Rajkiran
4.	09-05-2026	-	16	-	3	19	11-05-2026	10:10am	Saravanan	Rajkiran
5.	12-05-2026	-	18	-	-	18	13-05-2026	10.15am	Saravanan	Rajkiran
6.	14-05-2026	-	48	-	-	48	15-05-2026	10:30am	Saravanan	Rajkiran
7.	16-05-2026	-	21	-	-	21	18-05-2026	11:00am	Saravanan	Rajkiran
8.	19-05-2026	-	16	-	-	16	20-05-2026	11:00am	Saravanan	Rajkiran
9.	21-05-2026	-	48	-	-	48	22-05-2026	10:35am	Saravanan	Rajkiran
10.	23-05-2026	-	16	-	5	21	25-05-2026	11:00am	Saravanan	Rajkiran
11.	26-05-2026	-	42	-	-	42	27-05-2026	11:00am	Saravanan	Rajkiran
12.	28-05-2026	1	18	-	-	19	29-05-2026	10:35am	Saravanan	Rajkiran
13.										
14.										
	TOTAL (Kgs)	1	295	-	10	306				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF...

(Sign & Date)

Checked by

01 June 2026 | 4:44 AM EDT

DocuSigned by:

G. Hubert Vino

02CF50F7F1A44D8...

(Sign & Date)

Verified by

01 June 2026 | 5:39 AM EDT