



Biomedical Waste Disposal Record

RHS_F_GN_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: June 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.	31-05-2025	-	31	-	3	34	02-06-2025	10:21am	Saravanan	Deepak
2.	03-06-2025	-	18	-	-	18	04-06-2025	10:39am	Saravanan	Rajkiran
3.	05-06-2025	-	31	-	-	31	06-06-2025	10:59am	Saravanan	Rajkiran
4.	07-06-2025	-	27	-	-	27	09-06-2025	17:05pm	Magesh	Jintu
5.	10-06-2025	-	18	-	4	22	11-06-2025	11.00am	Saravanan	Rajkiran
6.	12-06-2025	-	33	-	1	34	13-06-2025	10:23am	Saravanan	Rajkiran
7.	16-06-2025	-	16	-	-	16	16-06-2025	10:50am	Saravanan	Jintu
8.	18-06-2025	-	54	-	-	54	18-06-2025	10:50am	Saravanan	Jintu
9.	20-06-2025	1	17	-	-	18	20-06-2025	10:35am	Saravanan	Jintu
10.	21-06-2025	-	18	-	5	23	23-06-2025	10:38am	Saravanan	Rajkiran
11.	24-06-2025	-	31	-	2	33	25-06-2025	11:05am	Saravanan	Rajkiran
12.	26-06-2025	-	27	-	-	27	27-06-2025	10:30am	Saravanan	Rajkiran
13.	30-06-2025	-	08	-	2	10	30-06-2025	11:10am	Saravanan	Jintu
14.										
	TOTAL (Kgs)	1	329	-	17	347				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF...

(Sign & Date)

Checked by

01 July 2025 | 4:41 AM EDT

DocuSigned by:

G. Hubert Vino

02CF50F7F1A44D8...

(Sign & Date)

Verified by

01 July 2025 | 4:45 AM EDT