



Biomedical Waste Disposal Record

RHS_F_GN_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: September 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.	30-08-2025	-	21	-	-	21	01-09-2025	10:35am	Saravanan	Rajkiran
2.	02-09-2025	-	18	-	-	18	03-09-2025	10:40am	Saravanan	Rajkiran
3.	04-09-2025	-	28	-	3	31	05-09-2025	11:15am	Saravanan	Rajkiran
4.	06-09-2025	-	37	-	2	39	08-09-2025	11:10pm	Saravanan	Jintu Das
5.	09-09-2025	-	47	-	-	47	10-09-2025	11.05am	Saravanan	Jintu Das
6.	11-09-2025	-	18	-	-	18	12-09-2025	10:55am	Saravanan	Jintu Das
7.	13-09-2025	-	27	-	2	29	15-09-2025	10:35am	Saravanan	Rajkiran
8.	16-09-2025	-	31	-	-	31	17-09-2025	11:05am	Saravanan	Rajkiran
9.	18-09-2025	-	18	-	4	22	19-09-2025	10:40am	Saravanan	Rajkiran
10.	20-09-2025	-	18	-	4	22	22-09-2025	11:00am	Saravanan	Jintu Das
11.	23-09-2025	-	16	-	-	17	24-09-2025	11:05am	Saravanan	Jintu Das
12.	25-09-2025	-	28	-	-	28	26-09-2025	10:45am	Saravanan	Jintu Das
13.	27-09-2025	-	51	-	2	53	29-09-2025	10:38am	Saravanan	Rajkiran
14.										
	TOTAL (Kgs)	-	358	-	17	375				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF... (Sign & Date)

Checked by

01 October 2025 | 2:08 AM EDT

DocuSigned by:

G. Hubert Kino

02CF50F7F1A44D8... (Sign & Date)

Verified by

01 October 2025 | 4:54 AM EDT