



Biomedical Waste Disposal Record

RHS_F_GN_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: December 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.	29-11-2025	-	24	-	-	24	01-12-2025	10:52am	Saravanan	Jintu Das
2.	02-12-2025	-	27	-	5	32	03-12-2025	11:00am	Saravanan	Jintu Das
3.	04-12-2025	-	36	-	-	36	05-12-2025	10:55am	Saravanan	Jintu Das
4.	06-12-2025	-	38	-	5	43	08-12-2025	11:30am	Saravanan	Rajkiran
5.	09-12-2025	-	24	-	-	24	10-12-2025	10.45am	Saravanan	Rajkiran
6.	11-12-2025	-	27	-	11	38	12-12-2025	10:40am	Saravanan	Rajkiran
7.	13-12-2025	-	19	-	4	23	15-12-2025	11:35am	Saravanan	Jintu Das
8.	16-12-2025	-	26	-	-	26	17-12-2025	10:44am	Saravanan	Jintu Das
9.	18-12-2025	-	14	-	-	14	19-12-2025	10:38am	Saravanan	Jintu Das
10.	20-12-2025	-	31	-	-	31	22-12-2025	10:26am	Saravanan	Rajkiran
11.	23-12-2025	-	19	-	2	21	24-12-2025	11:00am	Saravanan	Rajkiran
12.	24-12-2025	-	6	-	-	6	26-12-2025	10:30am	Saravanan	Rajkiran
13.	27-12-2025	-	17	-	-	17	29-12-2025	10:40am	Saravanan	Jintu Das
14.	30-12-2025	-	31	-	-	31	31-12-2025	10:40am	Saravanan	Jintu Das
	TOTAL (Kgs)	-	339	-	27	366				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF...

(Sign & Date)

Checked by

02 January 2026 | 12:17 AM EST

DocuSigned by:

G. Hubert Vino

02CF50F7F1A44D8...

(Sign & Date)

Verified by

02 January 2026 | 2:21 AM EST