



Biomedical Waste Disposal Record

RHS_F_GN_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: August 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.	01-08-2026	-	31	-	5	36	03-08-2026	11:06am	Saravanan	Jintu Das
2.	04-08-2026	-	16	-	-	16	05-08-2026	11:14am	Saravanan	Jintu Das
3.	06-08-2026	-	17	-	-	17	07-08-2026	10:40am	Saravanan	Jintu Das
4.	08-08-2026	-	28	-	3	31	10-08-2026	11:12am	Saravanan	Jintu Das
5.	11-08-2026	-	32	-	-	32	12-08-2026	11:05am	Saravanan	Rajkiran
6.	13-08-2026	-	27	-	-	27	14-08-2026	10:40am	Saravanan	Rajkiran
7.	14-08-2026	-	20	-	5	25	17-08-2026	10:55am	Aravindh	Jintu Das
8.	18-08-2026	-	15	-	-	15	19-08-2026	11:10am	Ganesh	Jintu Das
9.	20-08-2026	-	32	-	-	32	21-08-2026	11:10am	Saravanan	Jintu Das
10.	22-08-2026	-	45	-	-	45	24-08-2026	10:00am	Aravindh	Rajkiran
11.	25-08-2026	-	30	-	-	30	26-08-2026	10:00am	Giri	Rajkiran
12.	27-08-2026	-	31	-	3	34	28-08-2026	10:30am	Saravanan	Rajkiran
13.	29-08-2026	-	21	-	-	21	31-08-2026	11:05am	Saravanan	Jintu Das
14.										
	TOTAL (Kgs)	-	345	-	16	361				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF...

(Sign & Date)

Checked by

02 September 2026 | 2:18 AM EDT

DocuSigned by:

G. Hubert Kuo

02CF50F7F1A44D8...

(Sign & Date)

Verified by

02 September 2026 | 2:25 AM EDT