



Biomedical Waste Disposal Record

RHS_F_GN_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: July 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.	02-07-2025	-	18	-	-	18	02-07-2025	11:50am	Saravanan	Jintu Das
2.	04-07-2025	-	17	-	-	17	04-07-2025	10:25am	Saravanan	Jintu Das
3.	05-07-2025	-	14	-	2	16	07-07-2025	10:20am	Saravanan	Rajkiran
4.	08-07-2025	-	19	-	4	23	09-07-2025	11:30pm	Saravanan	Rajkiran
5.	11-07-2025	-	29	-	-	29	11-07-2025	09:55am	Saravanan	Jintu Das
6.	14-07-2025	-	17	-	2	19	14-07-2025	10:25am	Saravanan	Jintu Das
7.	16-07-2025	-	14	-	-	14	16-07-2025	10:50am	Saravanan	Jintu Das
8.	18-07-2025	-	26	-	-	26	18-07-2025	10:35am	Saravanan	Jintu Das
9.	19-07-2025	1	32	-	-	32	21-07-2025	10:26am	Saravanan	Rajkiran
10.	22-07-2025	-	10	-	-	10	23-07-2025	09:57am	Abin	Rajkiran
11.	24-07-2025	-	15	-	-	15	25-07-2025	09:58am	Saravanan	Rajkiran
12.	28-07-2025	-	25	-	10	35	28-07-2025	10:25am	Saravanan	Jintu Das
13.	30-07-2025	2	34	-	-	36	30-07-2025	10:55am	Saravanan	Jintu Das
14.										
	TOTAL (Kgs)	2	270	-	18	290				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF...

(Sign & Date)

Checked by

02 August 2025 | 2:21 AM EDT

DocuSigned by:

G. Hubert Kiro

02CF50F7F1A44D8...

(Sign & Date)

Verified by

02 August 2025 | 3:39 AM EDT