

revvity

Biomedical Waste Disposal Record

RHS_F_GN_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: March 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.	28-02-2026	-	46	-	-	46	02-03-2026	11:00am	Saravanan	Rajkiran
2.	03-03-2026	-	52	-	-	52	04-03-2026	11:15am	Saravanan	Rajkiran
3.	05-03-2026	-	24	-	-	24	06-03-2026	11:15am	Saravanan	Rajkiran
4.	07-03-2026	-	43	-	3	46	09-03-2026	12:35am	Saravanan	Jintu Das
5.	10-03-2026	-	34	-	-	34	11-03-2026	10.40am	Saravanan	Jintu Das
6.	12-03-2026	-	24	-	-	24	13-03-2026	10:50am	Saravanan	Jintu Das
7.	14-03-2026	-	18	-	-	18	16-03-2026	11:00am	Saravanan	Rajkiran
8.	17-03-2026	-	17	-	-	17	18-03-2026	11:00am	Saravanan	Rajkiran
9.	19-03-2026	-	42	-	-	42	20-03-2026	11:00am	Saravanan	Rajkiran
10.	21-03-2026	-	11	-	5	16	23-03-2026	11:00am	Saravanan	Jintu Das
11.	24-03-2026	-	46	-	-	46	25-03-2026	11:10am	Saravanan	Jintu Das
12.	26-03-2026	-	28	-	-	28	27-03-2026	11:00am	Saravanan	Jintu Das
13.	28-03-2026	-	15	-	-	15	30-03-2026	10:55am	Saravanan	Jintu Das
14.										
	TOTAL (Kgs)	-	400	-	8	408				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF...

(Sign & Date)

Checked by

01 April 2026 | 2:45 AM EDT

DocuSigned by:

G. Hubert Kuo

02CF50F7F1A44D8...

(Sign & Date)

Verified by

01 April 2026 | 7:53 AM EDT