

revvity

Biomedical Waste Disposal Record

RHS\_F\_GN\_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: February 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.	31-01-2026	-	20	-	-	20	02-02-2026	10:15am	Ragu	Rajkiran
2.	03-02-2026	-	26	-	-	26	04-02-2026	11:03am	Saravanan	Rajkiran
3.	05-02-2026	-	27	-	-	27	06-02-2026	10:45am	Saravanan	Rajkiran
4.	07-02-2026	-	21	-	-	21	09-02-2026	11:10am	Saravanan	Jintu Das
5.	10-02-2026	-	17	-	2	19	11-02-2026	11.10am	Saravanan	Jintu Das
6.	12-02-2026	-	16	-	-	16	13-02-2026	10:55am	Saravanan	Jintu Das
7.	14-02-2026	-	15	-	-	15	16-02-2026	10:20am	Saravanan	Rajkiran
8.	17-02-2026	-	37	-	4	41	18-02-2026	11:15am	Saravanan	Rajkiran
9.	19-02-2026	-	16	-	-	16	20-02-2026	11:00am	Saravanan	Rajkiran
10.	21-02-2026	-	18	-	3	21	23-02-2026	10:50am	Saravanan	Jintu Das
11.	24-02-2026	-	57	-	-	57	25-02-2026	11:15am	Saravanan	Jintu Das
12.	26-02-2026	-	18	-	3	21	27-02-2026	11:15am	Saravanan	Jintu Das
13.										
14.										
	<b>TOTAL (Kgs)</b>	-	<b>288</b>	-	<b>12</b>	<b>300</b>				

DocuSigned by:

Baranidaran S

DC2050D1B1DE4DF...

(Sign &amp; Date)

Checked by 02 March 2026 | 2:27 AM EST

DocuSigned by:

G. Hubert Kino

02CF50F7F1A44D8...

(Sign &amp; Date)

Verified by 02 March 2026 | 4:16 AM EST