

revvity	<i>Biomedical Waste Disposal Record</i>	RHS_F_GN_028	v.1
Effective Date: 01/January/2024		Approval Date: 18/February/2019	

Month/ Year: January 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.	31-12-2025	-	3	-	4	7	01-12-2026	11:00am	Saravanan	Jintu Das
2.	03-01-2026	-	34	-	-	34	03-12-2026	10:30am	Saravanan	Rajkiran
3.	06-12-2026	-	31	-	-	31	05-12-2026	10:45am	Saravanan	Rajkiran
4.	08-12-2026	-	53	-	-	53	08-12-2026	11:08am	Saravanan	Rajkiran
5.	10-12-2026	-	17	-	4	21	10-12-2026	11.03am	Saravanan	Jintu Das
6.	13-12-2026	-	18	-	-	18	12-12-2026	09:40am	Saravanan	Jintu Das
7.	17-12-2026	-	17	-	-	17	15-12-2026	10:40am	Saravanan	Rajkiran
8.	20-12-2026	-	28	-	5	33	17-12-2026	11:00am	Saravanan	Rajkiran
9.	22-12-2026	-	38	-	3	41	19-12-2026	11:00am	Saravanan	Rajkiran
10.	27-12-2026	-	31	-	-	31	22-12-2026	11:10am	Saravanan	Jintu Das
11.	29-12-2026	-	48	-	4	52	24-12-2026	11:00am	Saravanan	Jintu Das
12.										
13.										
14.										
	TOTAL (Kgs)	-	318	-	20	338				

DocuSigned by:
Baranidaran S
DC2050D1B1DE4DF...

(Sign & Date)

Checked by

02 February 2026 | 2:19 AM EST

DocuSigned by:
G. Hubert Kuo
02CF50F7F1A44D8...

(Sign & Date)

Verified by

02 February 2026 | 6:16 AM EST