

revvity

Biomedical Waste Disposal Record

RHS_F_GN_028

v.1

Effective Date: 01/January/2024

Approval Date: 18/February/2019

Month/ Year: August 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.	31-07-2025	-	17	-	-	17	01-08-2025	10:15am	Saravanan	Jintu Das
2.	02-08-2025	2	10	-	-	12	04-08-2025	09:55am	Abin	Rajkiran
3.	05-08-2025	-	13	-	-	13	06-08-2025	09:40am	Abin	Rajkiran
4.	07-08-2025	-	15	-	5	20	08-08-2025	09:45pm	Abin	Rajkiran
5.	09-08-2025	-	53	-	4	57	11-08-2025	11.10am	Saravanan	Rajkiran
6.	12-08-2025	-	9	-	-	9	13-08-2025	10:33am	Saravanan	Jintu Das
7.	16-08-2025	-	24	-	3	27	18-08-2025	10:15am	Saravanan	Rajkiran
8.	19-08-2025	-	24	-	-	24	20-08-2025	11:05am	Saravanan	Rajkiran
9.	21-08-2025	-	24	-	-	24	22-08-2025	11:30am	Ssaravanan	Rajkiran
10.	24-08-2025	-	17	-	2	19	25-08-2025	10:30am	Saravanan	Jintu Das
11.	28-08-2025	-	28	-	3	31	29-08-2025	11:10am	Saravanan	Jintu Das
12.										
13.										
14.										
	TOTAL (Kgs)	2	234	-	17	253				

DocuSigned by:
Baranidaran S
DC2050D1B1DE4DF...

(Sign & Date)

Checked by

01 September 2025 | 2:46 AM EDT

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

(Sign & Date)

Verified by

01 September 2025 | 3:41 AM EDT