

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

Month/ Year: January 2024

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.	02-01-2024	12	49			61	03-01-2024	10:24am	Saravanan	Deepak
2.	04-01-2024	11	48			59	05-01-2024	10:20am	Saravanan	Deepak
3.	06-01-2024		15			15	08-01-2024	09:41am	Raghu	Suman Banik
4.	09-02-2024		30			30	10-01-2024	09.20am	Raghu	Suman Banik
5.	11-01-2024		45			45	12-01-2024	09.40am	Raghu	Suman Banik
6.	18-02-2024	12	61			73	19-01-2024	09:55am	Saravanan	Deepak
7.	20-02-2024	6	17			23	22-01-2024	09:50am	Saravanan	Suman Banik
8.	23-01-2024	6	27			33	24-01-2024	10:15am	Saravanan	Suman Banik
9.	27-01-2024	6	17			23	29-01-2024	10:25am	Saravanan	Deepak
10.	30-01-2024	7	29		4	40	31-01-2024	10:15am	Saravanan	Deepak
11.										
12.										
13.										
	TOTAL (Kgs)	60	338	0	4	402				

DocuSigned by:
Dayasagari V
D650AD69608E40E (Sign & Date)

Checked by QA 02 February 2024 | 5:48 PM IST

DocuSigned by:
Prabhu Selvaraj
A51ABF0C7A5F4D7... (Sign & Date)

Verified by QM 02 February 2024 | 5:50 PM IST