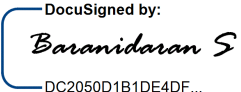
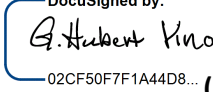


<b>revvity</b>	<i>Biomedical Waste Disposal Record</i>	<b>RHS_F_GN_028</b>	<b>v.1</b>
Effective Date: 01/January/2024		Approval Date: 18/February/2019	

**Month/ Year: March 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.	01-03-2025	-	16	-	1	17	03-03-2025	10:10am	Saravanan	Deepak
2.	04-03-2025	-	18	-	-	18	05-03-2025	10:20am	Saravanan	Deepak
3.	06-03-2025	-	49	-	-	49	07-03-2025	10:09am	Saravanan	Deepak
4.	08-03-2025	-	16	-	3	19	10-03-2025	10:40am	Saravanan	Deepak
5.	11-03-2025	-	21	-	-	21	12-03-2025	10:52am	Saravanan	Deepak
6.	13-03-2025	-	23	-	-	23	14-03-2025	09:52am	Abin	Deepak
7.	15-03-2025	-	32	-	-	32	17-03-2025	10:15am	Saravanan	Deepak
8.	18-03-2025	-	18	-	-	18	19-03-2025	10:27am	Saravanan	Deepak
9.	20-03-2025	-	24	-	1	25	21-03-2025	10:25am	Saravanan	Deepak
10.	22-03-2025	-	17	-	-	17	24-03-2025	10:20am	Saravanan	Deepak
11.	25-03-2025	-	18	-	4	22	26-03-2025	10:35am	Saravanan	Deepak
12.	27-03-2025	-	58	-	3	61	28-03-2025	10:30am	Saravanan	Deepak
13.	29-03-2025	-	15	-	-	15	31-03-2025	10:07am	Saravanan	Deepak
14.										
	<b>TOTAL (Kgs)</b>	-	<b>325</b>	-	<b>12</b>	<b>337</b>				

DocuSigned by:  
  
 DC2050D1B1DE4DF... (Sign & Date)  
**Checked by**  
 01 April 2025 | 5:54 AM EDT

DocuSigned by:  
  
 02CF50F7F1A44D8... (Sign & Date)  
**Verified by**  
 01 April 2025 | 10:26 AM EDT