
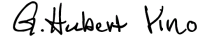


|                                 |   |                                 |            |
|---------------------------------|---|---------------------------------|------------|
| <b>revvity</b>                  | <i>Biomedical Waste Disposal Record</i> | <b>RHS_F_GN_028</b>             | <b>v.1</b> |
|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

**Month/ Year: October 2024**

| S. No. | Date of Generation | Quantity of BMW Generated (in KG)<br>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.     | 03-10-2024         | -   | 34         | -     | -        | 34         | 04-10-2024                   | 10:35am          | Saravanan               | Deepak        |
| 2.     | 05-10-2024         | -   | 31         | -     | 3        | 34         | 07-10-2024                   | 10:17am          | Saravanan               | Deepak        |
| 3.     | 08-10-2024         | -   | 31         | -     | -        | 31         | 09-10-2024                   | 10:40am          | Saravanan               | Deepak        |
| 4.     | 12-10-2024         | -   | 27         | -     | 2        | 29         | 14-10-2024                   | 10:50am          | Saravanan               | Deepak        |
| 5.     | 17-10-2024         | -   | 24         | -     | -        | 24         | 18-10-2024                   | 10.28am          | Saravanan               | Deepak        |
| 6.     | 19-10-2024         | -   | 26         | -     | -        | 26         | 21-10-2024                   | 11:30am          | Saravanan               | Deepak        |
| 7.     | 22-10-2024         | -   | 24         | -     | 2        | 26         | 23-10-2024                   | 10:35am          | Saravanan               | Deepak        |
| 8.     | 24-10-2024         | -   | 26         | -     | -        | 26         | 25-10-2024                   | 10:40am          | Saravanan               | Deepak        |
| 9.     | 26-10-2024         | -   | 14         | -     | -        | 14         | 28-10-2024                   | 10:05am          | Saravanan               | Deepak        |
| 10.    | 29-10-2024         | -   | 16         | -     | -        | 16         | 30-10-2024                   | 10:00am          | Saravanan               | Deepak        |
| 11.    |                    |   |            |       |          |            |                              |                  |                         |               |
| 12.    |                    |   |            |       |          |            |                              |                  |                         |               |
| 13.    |                    |   |            |       |          |            |                              |                  |                         |               |
| 14.    |                    |   |            |       |          |            |                              |                  |                         |               |
|        | <b>TOTAL (Kgs)</b> | -   | <b>253</b> | -     | <b>7</b> | <b>260</b> |                              |                  |                         |               |

DocuSigned by:  
  
 DC2050D1B1DE4DF... (Sign & Date)  
**Checked by**  
 01 November 2024 | 1:18 AM EDT

DocuSigned by:  
  
 02CF50F7F1A44D8... (Sign & Date)  
**Verified by**  
 01 November 2024 | 3:04 AM EDT