

**Resorufin, sodium salt \*CAS 34994-50-8\***Catalog number: 65  
Unit size: 100 mg**Product Details**

|                    |   |
|--------------------|---|
| Storage Conditions | Freeze (<-15 °C), Minimize light exposure |
| Expiration Date    | 24 months upon receiving                  |

**Chemical Properties**

|                    |   |
|--------------------|---|
| Appearance         | Dark purple solid   |
| Molecular Weight   | 235.17  |
| Soluble In         | Water   |
| Chemical Structure |  |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Excitation Wavelength | 571 nm |
| Emission Wavelength   | 584 nm |

**Applications**

Resorufin is used as a calibration standard for resorufin-derived enzyme substrates. It has pH dependent absorption and emission over the range of 5 to 9. It is also the reduction product of resazurin in cells.