

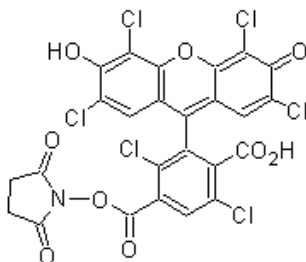
## Product Information Sheet

### Ordering Information

|                     |   |
|---------------------|---|
| Product Number:     | 202   |
| Product Name:       | 6-HEX, SE [6-Carboxy-2',4,4',5',7,7'-hexachlorofluorescein, succinimidyl ester] *Single Isomer* |
| Unit Size:          | 5 mg  |
| Storage Conditions: | <-15 °C and kept from light and moisture  |
| Expiration Date:    | 12 months upon receiving  |

### Chemical, Physical and Spectral Properties

|                     |           |
|---------------------|-----------|
| Molecular Weight:   | 680.06    |
| Appearance:         | Red solid |
| Chemical Structure: |           |



|                        |        |
|------------------------|--------|
| Soluble in:            | DMSO   |
| Excitation Wavelength: | 533 nm |
| Emission Wavelength:   | 550 nm |

### Application Notes

6-HEX, SE is the amine-reactive succinimidyl ester of 6-HEX acid. It is widely used in nucleic acid sequencing and related research.