

## 5(6)-Carboxyrhodamine 6G cadaverine

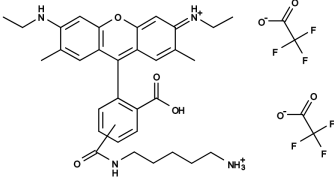
Catalog number: 343

Unit size: 1 mg

### Product Details

|                    |  |
|--------------------|--|
| Storage Conditions | Freeze (<math>< -15\text{ }^\circ\text{C}</math>), Minimize light exposure |
| Expiration Date    | 12 months upon receiving   |

### Chemical Properties

|                    |  |
|--------------------|--|
| Appearance         | Red solid  |
| Molecular Weight   | 770.72   |
| Soluble In         | DMSO   |
| Chemical Structure |  |

### Spectral Properties

|                       |        |
|-----------------------|--------|
| Excitation Wavelength | 522 nm |
| Emission Wavelength   | 546 nm |

### Applications

5(6)-Carboxyrhodamine 6G cadaverine is a building block for developing rhodamine bioconjugates. It is so a fluorescent transglutaminase substrate to label proteins by transamination. In general, rhodamine 6G conjugates have higher fluorescence quantum yields than tetramethylrhodamine conjugates, as well as good photostability. The excitation and emission wavelengths of rhodamine 6G fall between those of fluorescein and tetramethylrhodamine derivatives, making 5(6)-CR6G another color choice for the multicolor fluorescence imaging applications. 5(6)-CR6G is the mixture of two carboxy rhodamine 6G isomers.