

**Fluorescein-5-maleimide \*CAS 75350-46-8\***

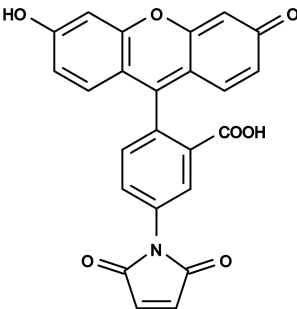
Catalog number: 130

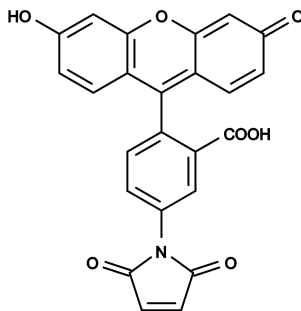
Unit size: 25 mg

**Product Details**

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

**Chemical Properties**

Appearance	Orange solid
Molecular Weight	427.36
Soluble In	DMSO
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	498 nm
Emission Wavelength	517 nm

**Applications**

Maleimides are among the most frequently used reagents for thiol modification. In most proteins, the site of reaction is at cysteine residues that either are intrinsically present or resulted from reduction of cystines. Unlike iodoacetamides, maleimides do not react with histidine and methionine under physiological conditions. Fluorescein-5-maleimide is one of the most popular fluorescent dyes for thiol modifications of proteins along with 5-TMRIA.