

# Acridinium Maleimide

Catalog Number: 26018

Unit Size: 1 mg

## Product Details

---

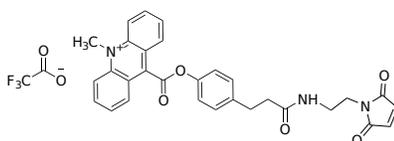
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

## Chemical Properties

---

Appearance	Solid
Molecular Weight	621.57
Soluble In	DMSO

Chemical Structure



## Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

## Applications

---

Acridinium is a univalent radical derived from acridine. Acridinium maleimides are synthetic molecules known for their strong chemiluminescence, commonly used for labeling thiol groups on proteins, oligonucleotide thiophosphates, or low molecular weight ligands. The chemiluminescence of acridinium maleimides is triggered by exposure to an alkaline hydrogen peroxide solution, facilitating their use in sensitive detection methodologies within immunoassays and various biological assays. For accurate signal detection in a chemiluminescence analyzer, a trigger solution containing (A) 0.1% H<sub>2</sub>O<sub>2</sub> and 0.1 M HNO<sub>3</sub>, and (B) 0.25 M NaOH and 2% Triton X-100 is recommended.