

# Norfentanyl Maleimide

Catalog Number: 50626

Unit Size: 1 mg

## Product Details

---

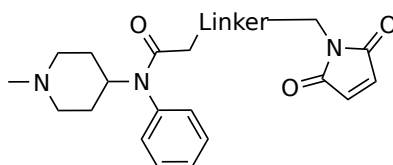
|                    |                                            |
|--------------------|--------------------------------------------|
| Storage Conditions | Freeze (< -15 °C), Minimize light exposure |
| Expiration Date    | 24 months upon receiving                   |

## Chemical Properties

---

|                  |        |
|------------------|--------|
| Appearance       | Solid  |
| Molecular Weight | 412.49 |
| Soluble In       | DMSO   |

Chemical Structure



## Spectral Properties

---

|                       |     |
|-----------------------|-----|
| Excitation Wavelength | N/A |
| Emission Wavelength   | N/A |

## Applications

---

Norfentanyl Maleimide is an excellent thiol-reactive building block for developing antibodies, assays and tests for detecting norfentanyl substances. It readily reacts with peptides, proteins and other thiol-containing biomolecules such as thiol-modified oligos. Norfentanyl is a monocarboxylic acid amide resulting from the formal condensation of the aryl amino group of 4-(N'-phenyl)piperidin-4-amine with propanoic acid. It is a major metabolite of fentanyl. It has a role as an opioid analgesic and a drug metabolite. It is an anilide, a member of piperidines and a monocarboxylic acid amide.