

# Norfentanyl-KLH Conjugate

Catalog Number: 50630

Unit Size: 1 mg

## Product Details

---

Storage Conditions Freeze (&lt; -15 °C), Minimize light exposure

Expiration Date 24 months upon receiving

## Chemical Properties

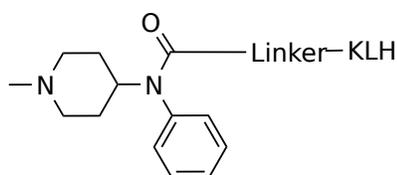
---

Appearance Solid

Molecular Weight N/A

Soluble In Water

Chemical Structure



## Spectral Properties

---

Excitation Wavelength N/A

Emission Wavelength N/A

## Applications

---

Norfentanyl-KLH Conjugate is an excellent antigen to develop anti-norfentanyl antibodies that can be used for detecting norfentanyl substances. 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.