

# Medetomidine azide

Catalog Number: 50652

Unit Size: 1 mg

## Product Details

---

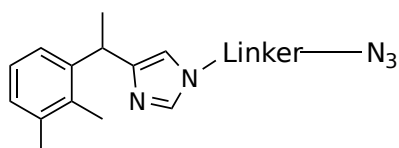
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

## Chemical Properties

---

Appearance	Solid
Molecular Weight	382.51
Soluble In	DMSO

Chemical Structure



## Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

## Applications

---

Medetomidine Azide is a functionalized derivative of medetomidine, a selective  $\alpha_2$ -adrenergic receptor agonist, modified to incorporate an azide (-N<sub>3</sub>) group for bioorthogonal conjugation via copper-catalyzed or copper-free click chemistry. The azide moiety allows for efficient and selective coupling to alkyne-functionalized molecules through Huisgen cycloaddition, enabling covalent attachment under mild conditions without interfering with biological activity. This reagent is ideal for applications requiring spatial and temporal control over conjugation, such as targeted drug delivery, receptor-ligand interaction studies, and the synthesis of functionalized probes or nanoparticles. Medetomidine Azide retains the pharmacological properties of the parent compound, making it a powerful tool for constructing receptor-targeted bioconjugates while preserving biological function.