

## EDTA Azide

Catalog Number: 65007

Unit Size: 5 mg

### Product Details

---

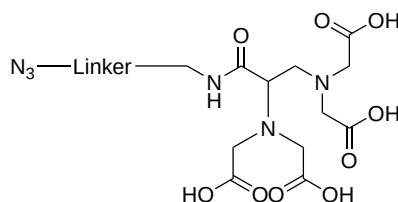
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

### Chemical Properties

---

Appearance	Solid colorless
Molecular Weight	768.77
Soluble In	DMF

Chemical Structure



### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

EDTA azide is an excellent building block to develop site-specific ethylenediaminetetraacetic acid (EDTA) probes for variety of biomedical applications. It readily reacts with an alkyne compound (such as DBCO) under the well-known click chemistry conditions. EDTA is the most commonly used chelator for complexing transitional metal ions. EDTA chelation therapy uses the chelating agent, EDTA, to remove heavy metals and minerals from the body. It involves injecting EDTA into the bloodstream, where it binds to these substances, allowing them to be excreted through urine. EDTA is approved for treating lead poisoning, but it's also used in other conditions, including coronary artery disease and arthritis.