

## 2',3'-cGAMP succinimidyl ester

Catalog Number: 20334

Unit Size: 100 ug

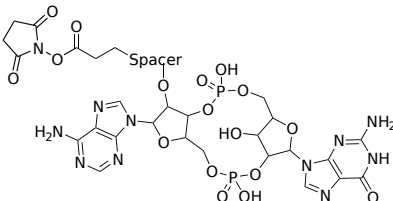
### Product Details

---

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

### Chemical Properties

---

Appearance	Solid colorless
Molecular Weight	1131.04
Soluble In	DMSO
Chemical Structure	

### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

2',3'-cGAMP succinimidyl ester is an excellent reagent to covalently conjugate 2',3'-cGAMP to biological targets, such as proteins, peptides, and nucleic acids. 2',3'-cGAMP has gained significant attention in recent years due to its potential as a therapeutic target for diseases such as cancer and viral infections. It has been shown to activate the immune system and enhance the efficacy of immune checkpoint inhibitors, which are a type of cancer immunotherapy. 2',3'-cGAMP (cyclic GMP-AMP) is a cyclic dinucleotide second messenger molecule that plays a critical role in the innate immune system. It is synthesized by the enzyme cGAS (cyclic GMP-AMP synthase) in response to cytosolic DNA that has been released from damaged or infected cells. Once synthesized, 2',3'-cGAMP binds to the adaptor protein STING (stimulator of interferon genes) and triggers downstream signaling pathways that result in the production of type I interferons and other cytokines, leading to an immune response against the invading pathogen.