

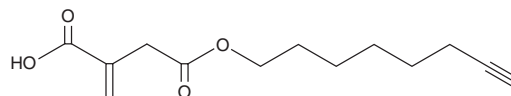
# PRODUCT INFORMATION



## 4-octynyl Itaconate

Item No. 28455

**CAS Registry No.:** 2454181-83-8  
**Formal Name:** 2-methylene-butanedioic acid,  
4-(7-octyn-1-yl) ester  
**Synonym:** Click Tag™ 4-octynyl Itaconate  
**MF:** C<sub>13</sub>H<sub>18</sub>O<sub>4</sub>  
**FW:** 238.3  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

4-octynyl Itaconate is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-octynyl itaconate in the solvent of choice, which should be purged with an inert gas. 4-octynyl Itaconate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 4-octynyl itaconate in these solvents is approximately 30 mg/ml.

### Description

4-octynyl Itaconate is an alkyne derivative of the itaconate prodrug 4-octyl itaconate (Item No. 25374). 4-octynyl Itaconate can be used for click chemistry applications.

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/25/2023

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM