

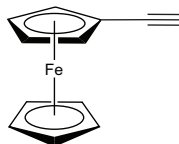
# PRODUCT INFORMATION



## Ethynylferrocene

Item No. 43793

CAS Registry No.: 1271-47-2  
Formal Name: ethynyl-ferrocene  
MF:  $C_{12}H_{10}Fe$   
FW: 210.1  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

Ethynylferrocene is supplied as a solid. A stock solution may be made by dissolving the ethynylferrocene in the solvent of choice, which should be purged with an inert gas. Ethynylferrocene is soluble ( $\geq 10$  mg/ml) in ethanol and DMSO.

### Description

Ethynylferrocene is a clickable form of the redox catalyst ferrocene.<sup>1</sup> It has been used to label ssDNA to measure DNA hybridization rates and to immobilize oligonucleotides to silicon-modified surfaces to detect complementary DNA strands.<sup>1,2</sup>

### References

1. Hu, Q., Deng, X., Kong, J., *et al.* Simple and fast electrochemical detection of sequence-specific DNA via click chemistry-mediated labeling of hairpin DNA probes with ethynylferrocene. *Analyst* 140(12), 4154-4161 (2015).
2. Ciampi, S., James, M., Michaels, P., *et al.* Tandem "click" reactions at acetylene-terminated Si(100) monolayers. *Langmuir* 27(11), 6940-6949 (2011).

#### 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, 06/24/2025

#### 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