

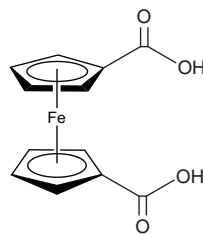
PRODUCT INFORMATION



1,1'-Ferrocenedicarboxylic Acid

Item No. 43795

CAS Registry No.: 1293-87-4
Formal Name: 1,1'-dicarboxy-ferrocene
Synonyms: DFc, $\text{Fc}(\text{COOH})_2$
MF: $\text{C}_{12}\text{H}_{10}\text{FeO}_4$
FW: 274.1
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

1,1'-Ferrocenedicarboxylic acid is supplied as a solid. A stock solution may be made by dissolving the 1,1'-ferrocenedicarboxylic acid in the solvent of choice, which should be purged with an inert gas. 1,1'-Ferrocenedicarboxylic acid is soluble (≥ 10 mg/ml) in DMSO.

Description

1,1'-Ferrocenedicarboxylic acid is a redox catalyst.^{1,2} It has been used in lithium ion-based dual-ion batteries.¹ 1,1'-Ferrocenedicarboxylic acid has also been conjugated to Prussian blue derivatives for the detection of hydrogen peroxide and glucose.²

References

1. Li, C., Yang, H., Xie, J., *et al.* Ferrocene-based mixed-valence metal-organic framework as an efficient and stable cathode for lithium-ion-based dual-ion battery. *ACS Appl. Mater. Interfaces* **12**(29), 32719-32725 (2020).
2. Tong, Y., Liu, Q., Yang, D., *et al.* Ferrocene derivatives constructed Prussian blue analogs for colorimetric detection of hydrogen peroxide and glucose. *Biosens. Bioelectron.* **287**, 117714 (2025).

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, 08/22/2025

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