

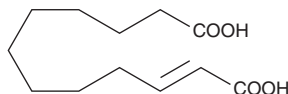
PRODUCT INFORMATION



Traumatic Acid

Item No. 88820

CAS Registry No.: 6402-36-4
Formal Name: 2E-dodecenedioic acid
Synonyms: FA 12:2;O2,
trans-2-Dodecenedioic acid
MF: C₁₂H₂₀O₄
FW: 228.3
Purity: ≥98%
Supplied as: A crystalline 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

Traumatic acid is supplied as a crystalline solid. Stock solutions of traumatic acid can be prepared by dissolving the traumatic acid in the solvent of choice, which should be purged with an inert gas. Traumatic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of traumatic acid in these solvents is approximately 10 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of traumatic acid is needed, it can be prepared by dissolving the crystalline compound in aqueous buffers. Traumatic acid is soluble in PBS (pH 7.2) at a concentration of approximately 900 µg/ml. The solubility of traumatic acid is increased in basic buffers. The solubility of traumatic acid in PBS (pH 9.0) is approximately 1.2 mg/ml. Store aqueous solutions of traumatic acid on ice and use within 12 hours of preparation. Although the aqueous solutions of traumatic acid may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

Description

Traumatic acid is a product of the hydroperoxide lyase pathway in plants. Traumatic acid is a wound-healing agent that stimulates cell division near a wound site to form a protective callus.¹

Reference

1. Vick, B.A. *Oxygenated fatty acids of the lipoxygenase pathway*. In: Lipid Metabolism in Plants (Moore Jr., T.S., ed.) pp. 167-191, CRC Press, Boca Raton (1993).

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, 02/09/2024

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