

PRODUCT INFORMATION



DL-erythro/threo-Isocitric Acid (sodium salt hydrate)

Item No. 32541

Formal Name: 1-hydroxypropane-1,2,3-tricarboxylate, trisodium salt hydrate

MF: $C_6H_5O_7 \cdot 3Na [XH_2O]$

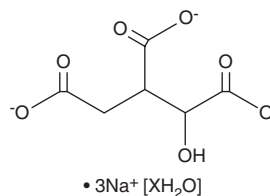
FW: 258.1

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years



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

Laboratory Procedures

DL-erythro/threo-Isocitric acid (sodium salt hydrate) is supplied as a solid. Aqueous solutions of DL-erythro/threo-isocitric acid can be prepared by directly dissolving the DL-erythro/threo-isocitric acid in aqueous buffers. The solubility of DL-erythro/threo-isocitric acid in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

DL-erythro/threo-Isocitric acid is a mixture of D-erythro-isocitric acid, L-erythro-isocitric acid, D-threo-isocitric acid, and L-threo-isocitric acid.

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, 09/29/2022

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