

Product Information



Aconitase Isocitric Dehydrogenase

Catalog No. 705510

Laboratory Procedures

The vial contains lyophilized isocitric dehydrogenase. Aconitase Isocitric Dehydrogenase, when stored as supplied, is stable for six months at -20°C. Prior to use in the Aconitase Assay Kit (Catalog No. 705502), reconstitute the contents of the vial with 2 ml of HPLC-grade water and store on ice. The reconstituted isocitric dehydrogenase is stable for one hour on ice. One vial of isocitric dehydrogenase is sufficient to evaluate 32 wells.

CAUTIONS

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work exclusively with the Aconitase Assay Kit from Cayman Chemical (Catalog No. 705502). This product may not perform as described if used with other assay reagents or protocols.

Related Products

96-Well Cover Sheets - Cat. No. 400012 • 96-Well Plate (Colorimetric Assay) - Cat. No. 400014 • Aconitase Assay Kit - Cat. No. 705502 • Aconitase Assay Buffer (10X) - Cat. No. 705503 • Aconitase Positive Control - Cat. No. 705504 • Aconitase Inhibitor - Cat. No. 705505 • Aconitase Substrate Solution (10X) - Cat. No. 705506 • Aconitase Cysteine Hydrochloride - Cat. No. 705507 • Aconitase Ferrous Ammonium Sulfate - Cat. No. 705508 • Aconitase NADP⁺ Reagent - Cat. No. 705509

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSD supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 10/21/2008