

PRODUCT INFORMATION



Krebs Cycle Standard Mixture

Item No. 28681

Supplied as: A solution in water

Storage: 4°C

Stability: ≥1 year

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

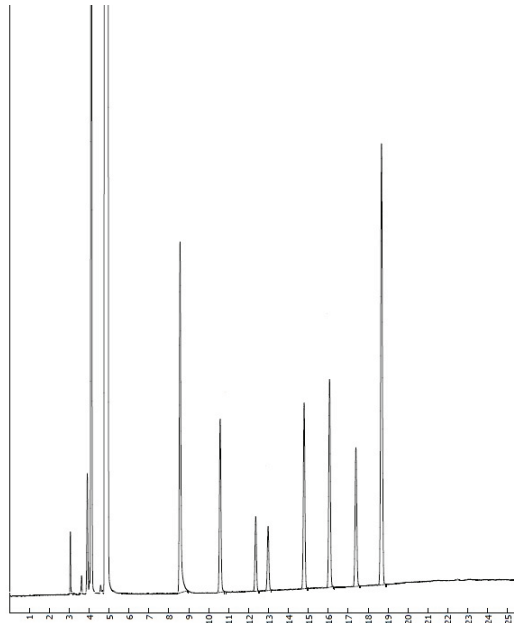
Description and Data

Krebs Cycle Standard Mixture contains seven non-volatile Krebs cycle metabolic intermediates. It is intended for use as a reference standard for the detection of these intermediates in various sample types by GC-FID, GC-MS, or LC-MS. For some applications, derivatization may be required.

Included are:

Elution Order*	Item Name
1	Pyruvic Acid
2	Lactic Acid
3	Oxaloacetic Acid
4	Oxalic Acid
5	Methyl Malonic Acid
6	Fumaric Acid
7	Succinic Acid

*GC Column Used: Supelcowax, 30 m x 0.53 mm x 0.2 µm



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, 03/31/2026

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