



CERTIFICATE OF ACCREDITATION

The ANSI National Accreditation Board

Hereby attests that

Cayman Chemical Company Incorporated

1180 East Ellsworth Road
Ann Arbor, MI 48108

Fulfills the requirements of

ISO/IEC 17025:2017

In the field of

TESTING

This certificate is valid only when accompanied by a current scope of accreditation document.
The current scope of accreditation can be verified at www.anab.org.

Jason Stine, Vice President

Expiry Date: 21 April 2027

Certificate Number: AT-1773



This laboratory is accredited in accordance with the recognized International Standard ISO/IEC 17025:2017.
This accreditation demonstrates technical competence for a defined scope and the operation of a laboratory
quality management system (refer to joint ISO-ILAC-IAF Communiqué dated April 2017).

SCOPE OF ACCREDITATION TO ISO/IEC 17025:2017

Cayman Chemical Company Incorporated

1180 East Ellsworth Road

Ann Arbor, MI 48108

Roxanne Franckowski (734)975-3975

rfranckowski@caymanchem.com www.caymanchem.com

TESTING

Valid to: **April 21, 2027**

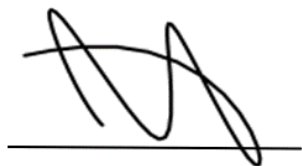
Certificate Number: **AT-1773**

Chemical

Specific Tests and/or Properties Measured	Specification, Standard, Method, or Test Technique	Items, Materials or Product Tested	Key Equipment or Technology
Reference Materials and Certified Reference Materials-Identifications and/or Concentrations	Internal SOP's Reference Methods: USP <621>, <736>, <854>	Reference Materials and Certified Reference Materials Organic Chemicals Neat and Dilute Materials Single and multi-component organic materials in solution or matrix	• GC/FID • GC/MS • HPLC-DAD • HPLC-MS • Residue on Ignition • FTIR • Loss on Drying • Water Determination (Karl Fischer Analysis) • Residual Solvent • Headspace Analysis • Thermogravimetric Analysis • qNMR

Note:

1. This scope is formatted as part of a single document including Certificate of Accreditation No. AT-1773.



Jason Stine, Vice President