



# CERTIFICATE OF ACCREDITATION

## ANSI-ASQ National Accreditation Board

500 Montgomery Street, Suite 625, Alexandria, VA 22314, 877-344-3044

This is to certify that

**Cayman Chemical Company**  
**1180 East Ellsworth Rd**  
**Ann Arbor, MI 48108**

has been assessed by ANAB  
and meets the requirements of international standard

## ISO/IEC 17025:2005

while demonstrating technical competence in the fields of

## TESTING

Refer to the accompanying Scope of Accreditation for information regarding the types of calibrations and/or tests to which this accreditation applies.

AT-1773

Certificate Number

  
ANAB Approval

Certificate Valid: 03/03/2018-04/21/2019  
Version No. 005 Issued: 03/03/2018



This laboratory is accredited in accordance with the recognized International Standard ISO/IEC 17025:2005. 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 17025:2005**

**Cayman Chemical Company**

1180 East Ellsworth Rd, Ann Arbor, MI 48108

Roxanne Franckowski 734-975-3975

[rfranckowski@caymanchem.com](mailto:rfranckowski@caymanchem.com) [www.caymanchem.com](http://www.caymanchem.com)

**TESTING**

Valid to: **April 21, 2019**

Certificate Number: **AT-1773**

**Organic Chemistry**

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>, <741>, <851>	Reference Materials and Certified Reference Materials Organic Chemicals Neat and Dilute Materials Single and multi-component organic materials in solution or matrix	<ul style="list-style-type: none"> <li>• GC/FID</li> <li>• GC/MS</li> <li>• HPLC-DAD</li> <li>• HPLC-MS</li> <li>• Residue on Ignition</li> <li>• FTIR</li> <li>• Loss on Drying</li> <li>• Water Determination (Karl Fischer Analysis)</li> <li>• Melting Point</li> <li>• Residual Solvent Headspace Analysis</li> </ul>

Note:

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



\_\_\_\_\_  
Vice President