



# CERTIFICATE OF ACCREDITATION

## ANSI National Accreditation Board

11617 Coldwater Road, Fort Wayne, IN 46845 USA

This is to certify that

### Cayman Chemical Company

1180 East Ellsworth Road

Ann Arbor, MI 48108

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

## ISO/IEC 17025:2017

while demonstrating technical competence in the field of

## TESTING

Refer to the accompanying Scope of Accreditation for information regarding the types of activities to which this accreditation applies

AT-1773

Certificate Number

  
ANAB Approval

Certificate Valid Through: 04/21/2021  
Version No. 007      Issued: 03/06/2020



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

1180 East Ellsworth Road

Ann Arbor, MI 48108

Roxanne Franckowski (734)975-3975

rfranckowski@caymanchem.com www.caymanchem.com

### TESTING

Valid to: **April 21, 2021**

Certificate Number: **AT-1773**

#### 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>, <854>	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>• Residual Solvent</li> <li>• Headspace Analysis</li> <li>• Thermogravimetric Analysis</li> </ul>

Note:

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



Vice President