



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. 006 Issued: 02/27/2019



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"> GC/FID GC/MS HPLC-DAD HPLC-MS Residue on Ignition FTIR Loss on Drying Water Determination (Karl Fischer Analysis) Melting Point Residual Solvent Headspace Analysis

Note:

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



 Vice President

