



# 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 17034:2016

while demonstrating technical competence in the field of

## REFERENCE MATERIAL PRODUCER

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

AR-1774

Certificate Number



ANAB Approval

Certificate Valid Through: 04/21/2021  
Version No. 007 Issued: 02/27/2019





SCOPE OF ACCREDITATION TO ISO 17034:2016

Cayman Chemical Company

1180 East Ellsworth Road

Ann Arbor, MI 48108

Roxanne Franckowski (734) 975-3975

rfranckowski@caymanchem.com www.caymanchem.com

REFERENCE MATERIAL PRODUCER

Valid to: April 21, 2021

Certificate Number: AR-1774

Chemistry

Sub-Category of Reference Material	ILAC RM Category	Class or Type of Reference Materials Produced (Include Range Where Applicable)	Methods or Techniques Used in the RMP Laboratory (if Appropriate)
Reference Materials and Certified Reference Materials	A 3.1	Single component or multi-component organic materials either neat or in dilute organic or aqueous solvents.	<ul style="list-style-type: none"> <li>GC/FID, USP &lt;621&gt;</li> <li>GC/MS, USP &lt;736&gt;</li> <li>HPLC-DAD, USP &lt;621&gt;</li> <li>HPLC-MS, USP &lt;736&gt;</li> <li>Residue on Ignition, Cayman method ISO TST SD24</li> </ul>
Forensic Reference Materials and Certified Reference Materials	A 8	CRM Categories: <ul style="list-style-type: none"> <li>Drugs of Abuse</li> <li>Metabolites</li> <li>Stable Isotope labeled materials</li> <li>Impurities and Degradants</li> <li>Intermediates</li> <li>Pharmaceuticals</li> <li>Phytochemicals</li> </ul>	<ul style="list-style-type: none"> <li>FTIR, USP &lt;854&gt;</li> <li>Loss on Drying, Cayman method ISO TST SD06</li> <li>Water Determination (Karl Fischer Analysis), Cayman method ISO TST SD02</li> <li>Melting Point, USP &lt;741&gt;</li> <li>Residual Solvent Headspace Analysis, Cayman method ISO TST SD11</li> </ul>
Pure Organic Compounds			

Notes:

- Please contact the RMP organization for more information on CRM uncertainty values, Ucrm values, and other specific lot values. Some of this information may also be available on the RMP's website.
- This scope is formatted as part of a single document including Certificate of Accreditation No. AR-1774.



Vice President

