



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. 008 Issued: 03/06/2020



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"> GC/FID, USP <621> GC/MS, USP <736> HPLC-DAD, USP <621> HPLC-MS, USP <736> Residue on Ignition, Internal SOP FTIR, USP <854> Loss on Drying, Internal SOP Water Determination (Karl Fischer Analysis), Internal SOP Residual Solvent Headspace Analysis, Internal SOP Thermogravimetric Analysis, Internal SOP
Forensic Reference Materials and Certified Reference Materials	A 8	CRM Categories: <ul style="list-style-type: none"> Drugs of Abuse Metabolites Stable Isotope labeled materials Impurities and Degradants Intermediates Pharmaceuticals Phytochemicals 	
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