



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 Guide 34:2009

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 03/03/2018-04/21/2019
Version No. 006 Issued: 03/03/2018



SCOPE OF ACCREDITATION TO ISO GUIDE 34:2009

Cayman Chemical Company

1180 East Ellsworth Rd

Ann Arbor, MI 48108

Roxanne Franckowski 734-975-3975

rfranckowski@caymanchem.com www.caymanchem.com

REFERENCE MATERIAL PRODUCER

Valid to: **April 21, 2019**

Certificate Number: **AR-1774**

I. Chemical

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 2.7 A 8.2 A 2.6	Single component or multi-component organic materials either neat or in dilute organic or aqueous solvents. CRM Categories: <ul style="list-style-type: none"> • Drugs of Abuse • Metabolites • Stable Isotope labeled materials • Impurities and Degradants • Intermediates • Pharmaceuticals 	<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

Notes:

1. 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.
2. This scope is formatted as part of a single document including Certificate of Accreditation No. AR-1774.



Vice President