



# CERTIFICATE OF ACCREDITATION

## The ANSI National Accreditation Board

Hereby attests that

**Cayman Chemical Company Incorporated**  
**1180 East Ellsworth Road**  
**Ann Arbor, MI 48108**

Fulfills the requirements of

**ISO 17034:2016**

In the field of

**REFERENCE MATERIAL PRODUCER**

This certificate is valid only when accompanied by a current scope of accreditation document.  
The current scope of accreditation can be verified at [www.anab.org](http://www.anab.org).

A handwritten signature in black ink, appearing to be 'Jason Stine', is positioned above a horizontal line.

Jason Stine, Vice President

Expiry Date: 21 April 2027

Certificate Number: AR-1774



This reference material producer is accredited in accordance with the recognized International Standard ISO 17034:2016.  
This accreditation demonstrates technical competence for a defined scope and the operation of a reference material producer quality management system.

**SCOPE OF ACCREDITATION TO ISO 17034:2016**

**Cayman Chemical Company Incorporated**

1180 East Ellsworth Road

Ann Arbor, MI 48108

Roxanne Franckowski (734) 975-3975

[rfranckowski@caymanchem.com](mailto:rfranckowski@caymanchem.com) [www.caymanchem.com](http://www.caymanchem.com)

**REFERENCE MATERIAL PRODUCER**

Valid to: **April 21, 2027**

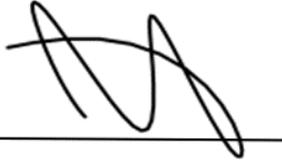
Certificate Number: **AR-1774**

**Chemical Properties**

Type of Reference Material	Description of the Reference Material Matrix or Artifact including the Property-Properties Characterized	Method or Techniques Used by the RMP Laboratory to Determine the Assigned Value (if Appropriate)
Reference Materials and Certified Reference Materials	<p style="text-align: center;">Forensic &amp; Pure Organic Compounds</p> <p>Single component or multi-component organic materials either neat or in dilute organic or aqueous solvents.</p> <p>CRM Categories:</p> <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>	<p>Characterization:</p> <ul style="list-style-type: none"> <li>• Characterization of non-operationally defined measurand using two or more methods of demonstrable accuracy in one or more competent laboratories</li> <li>• Value transfer from an RM to a closely matched candidate RM performed using a single measurement procedure performed by one laboratory</li> <li>• Characterization based on mass or volume of ingredients used in the preparation of the RM</li> <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 Headspace Analysis</li> <li>• Thermogravimetric Analysis</li> <li>• qNMR</li> </ul>

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.



Jason Stine, Vice President

