

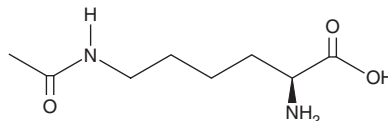
# PRODUCT INFORMATION



## L-Lys(Ac)-OH

Item No. 44219

CAS Registry No.: 692-04-6  
Formal Name: N<sup>6</sup>-acetyl-L-lysine  
Synonyms: AcK, N-ε-acetyl-L-Lysine  
MF: C<sub>8</sub>H<sub>16</sub>N<sub>2</sub>O<sub>3</sub>  
FW: 188.2  
Purity: ≥98%  
Supplied as: A solid  
Storage: -20°C  
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

L-Lys(Ac)-OH is supplied as a solid. Aqueous solutions of L-Lys(Ac)-OH can be prepared by directly dissolving the solid in aqueous buffers. L-Lys(Ac)-OH is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

L-Lys(Ac)-OH is an acetylated form of L-lysine (Item No. 36455). It is increased in the plasma of patients with primary biliary cholangitis- or primary sclerosing cholangitis-associated liver cirrhosis and is correlated with disease severity.<sup>1</sup>

### Reference

1. Mindikoglu, A.L., Coarfa, C., Opekun, A.R., *et al.* Metabolomic biomarkers are associated with mortality in patients with cirrhosis caused by primary biliary cholangitis or primary sclerosing cholangitis. *Future Sci. OA.* **6**(2), FSO441 (2019).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/03/2025

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM