

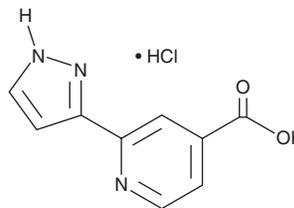
# PRODUCT INFORMATION



## 2-(1H-Pyrazol-3-yl)pyridine-4-carboxylic Acid (hydrochloride)

Item No. 43958

**CAS Registry No.:** 2749892-97-3  
**Formal Name:** 2-(1H-pyrazol-3-yl)-4-pyridinecarboxylic acid, monohydrochloride  
**MF:** C<sub>9</sub>H<sub>7</sub>N<sub>3</sub>O<sub>2</sub> • HCl  
**FW:** 225.6  
**Purity:** ≥95%  
**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

2-(1H-Pyrazol-3-yl)pyridine-4-carboxylic acid (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the 2-(1H-pyrazol-3-yl)pyridine-4-carboxylic acid (hydrochloride) in the solvent of choice, which should be purged with an inert gas. 2-(1H-Pyrazol-3-yl)pyridine-4-carboxylic acid (hydrochloride) is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

### Description

2-(1H-Pyrazol-3-yl)pyridine-4-carboxylic acid is a heterocycle that has been used as a cofactor in crystallography studies of lysine-specific demethylase 4A (KDM4A), also known as Jumanji domain containing 2A (JMJD2A).<sup>1</sup>

### Reference

1. Giri, N.C., Passantino, L., Sun, H., *et al.* Structural investigations of the nickel-induced inhibition of truncated constructs of the JMJD2 family of histone demethylases using X-ray absorption spectroscopy. *Biochemistry* **52**(24), 4168-4183 (2013).

#### 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, 08/11/2025

#### 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