

---

## 2,4 pyridindicarboxylic acid, hydrate

(Catalog No. M55002)

---

### Alternative Names

Lutidinic acid, PD2, AC1Q5UFW, pyridine carboxylate

### Description

Inhibitor of 2-OG-dependent histone demethylase JMJD2E, and trimethyllysine-specific histone demethylase JMJD2A.

### CAS Number

207671-42-9

### Chemical Formula

$C_7H_5NO_4 \cdot H_2O$

### Molecular Weight

167.1

### Size

1 g

### Purity

>98%

### Solubility

Water

### Storage Temp

-20°C

### Ordering Information:

Products	Size	Cat. No.
2,4 pyridindicarboxylic acid, hydrate	1 g	M55002 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*