

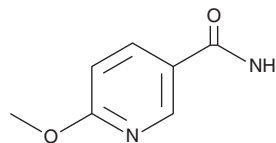
# PRODUCT INFORMATION



**JBSNF-000088**

Item No. 29920

**CAS Registry No.:** 7150-23-4  
**Formal Name:** 6-methoxy-3-pyridinecarboxamide  
**Synonyms:** 6-Methoxynicotinamide, NSC 70628  
**MF:** C<sub>7</sub>H<sub>8</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 152.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 240, 273 nm  
**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

JBSNF-000088 is supplied as a solid. A stock solution may be made by dissolving the JBSNF-000088 in the solvent of choice, which should be purged with an inert gas. JBSNF-000088 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JBSNF-000088 in ethanol is approximately 1 mg/ml and approximately 5 mg/ml in DMSO and DMF.

JBSNF-000088 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JBSNF-000088 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. JBSNF-000088 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

JBSNF-000088 is an inhibitor of nicotinamide N-methyltransferase (NNMT; IC<sub>50</sub>s = 1.8, 2.8, and 5 μM for human, monkey, and mouse NNMT, respectively).<sup>1</sup> It inhibits NNMT and reduces 1-methyl-nicotinamide (MNA) levels in U2OS and 3T3L1 cells (IC<sub>50</sub>s = 1.6 and 6.3 μM, respectively). JBSNF-000088 (50 mg/kg) reduces visceral white adipose tissue (WAT) MNA levels, body weight, fed blood glucose levels, and plasma and liver triglyceride levels, and improves oral glucose tolerance in a mouse model of diet-induced obesity (DIO). It also improves glucose tolerance, without affecting body weight, in the *ob/ob* and *db/db* mouse models of insulin resistance and diabetes, respectively.

## Reference

1. Kannt, A., Rajagopal, S., Kadnur, S.K., *et al.* A small molecule inhibitor of nicotinamide N-methyltransferase for the treatment of metabolic disorders. *Sci. Rep.* **8**(1), 3660 (2018).

### 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, 11/28/2022

## 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](http://WWW.CAYMANCHEM.COM)