

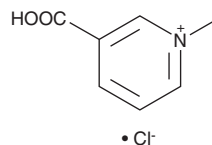
PRODUCT INFORMATION



Trigonelline (chloride)

Item No. 11904

CAS Registry No.: 6138-41-6
Formal Name: 3-carboxy-1-methyl-pyridinium, monochloride
Synonym: N-methyl Nicotinic Acid Betaine
MF: $C_7H_8NO_2 \cdot Cl$
FW: 173.6
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 265 nm
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years
Item Origin: Plant/*Trigonella foenum*



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

Laboratory Procedures

Trigonelline (chloride) is supplied as a crystalline solid. Aqueous solutions of trigonelline (chloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of trigonelline (chloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Trigonelline is an alkaloid that has been found in *L. japonicus* and an active metabolite of niacin that has diverse biological activities.¹⁻⁴ It increases levels of the brown fat marker proteins PRDM16, PGC-1 α , and UCP1 in 3T3-L1 adipocytes when used at a concentration of 75 μM .¹ Trigonelline inhibits degranulation of, and decreases the production of IL-6 and TNF- α in, activated primary mouse bone marrow mast cells (BMMCs).² *In vivo*, trigonelline (200 mg/kg) reduces serum IgE levels, pulmonary immune cell infiltration, and mucus secretion in a mouse model of ovalbumin-induced allergic asthma. It reduces serum levels of IL-1 β , IL-6, IL-18, and malondialdehyde (MDA) and renal cell apoptosis, as well as increases protein levels of peroxisome proliferator-activated receptor γ (PPAR γ) in a rat model of high-fat diet- and streptozotocin-induced type 2 diabetic nephropathy when administered at a dose of 40 mg/kg.³ Trigonelline (50 mg/kg) reduces hepatic *de novo* lipogenesis, induces hepatic autophagy, and prevents weight gain, insulin resistance, and hepatic steatosis in a mouse model of high-cholesterol and high-fat diet-induced non-alcoholic fatty liver disease (NAFLD).⁴

References

1. Choi, M., Mukherjee, S., and Yun, J.W. Trigonelline induces browning in 3T3-L1 white adipocytes. *Phytother Res.* **35**(2), 1113-1124 (2021).
2. Zhang, W., Zhang, Y., Chen, S., et al. Trigonelline, an alkaloid from *Leonurus japonicus* Houtt., suppresses mast cell activation and OVA-induced allergic asthma. *Front. Pharmacol.* **12**, 687970 (2021).
3. Li, Y., Li, Q., Wang, C., et al. Trigonelline reduced diabetic nephropathy and insulin resistance in type 2 diabetic rats through peroxisome proliferator-activated receptor- γ . *Exp. Ther. Med.* **18**(2), 1331-1337 (2019).
4. Sharma, L., Lone, N.A., Knott, R.M., et al. Trigonelline prevents high cholesterol and high fat diet induced hepatic lipid accumulation and lipo-toxicity in C57BL/6J mice, via restoration of hepatic autophagy. *Food Chem. Toxicol.* **121**, 283-296 (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, 01/04/2023

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