

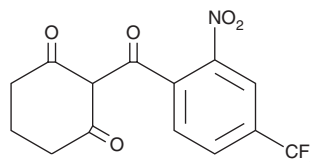
PRODUCT INFORMATION



Nitisinone

Item No. 17924

CAS Registry No.: 104206-65-7
Formal Name: 2-[2-nitro-4-(trifluoromethyl)benzoyl]-1,3-cyclohexanedione
Synonyms: NTBC, SC-0735
MF: C₁₄H₁₀F₃NO₅
FW: 329.2
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 271 nm
Supplied as: A crystalline 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

Nitisinone is supplied as a crystalline solid. A stock solution may be made by dissolving the nitisinone in the solvent of choice, which should be purged with an inert gas. Nitisinone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of nitisinone in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Nitisinone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nitisinone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Nitisinone has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Nitisinone is an inhibitor of 4-hydroxyphenylpyruvate dioxygenase (HPPD), which converts 4-hydroxyphenylpyruvate (HPPA) to homogentisate in the tyrosine catabolic pathway.¹ Nitisinone increases urinary levels of HPPA and 4-hydroxyphenyllactate (HPLA) in rats when administered at a dose of 10 mg/kg. Nitisinone (3 mg/kg) prevents the neonatal lethality of fumarylacetoacetate hydrolase (FAH) deficiency in mice when administered to pregnant dams.² It exhibits hepatoprotective effects in FAH^{-/-} mice, such as prevention of increases in plasma levels of aspartate serine aminotransferase (AST) and conjugated bilirubin, when administration is continued following birth at a dose of 1 mg/kg. Nitisinone (100 μg) decreases urinary excretion of homogentisate and increases urinary excretion of HPPA, HPLA, and 4-hydroxyphenylacetate in a mouse model of alkaptonuria induced by ethylnitrosourea.³ Formulations containing nitisinone have been used in the treatment of hereditary tyrosinemia type 1 (HT-1).

References

1. Ellis, M.K., Whitfield, A.C., Gowans, L.A., *et al.* Inhibition of 4-hydroxyphenylpyruvate dioxygenase by 2-(2-nitro-4-trifluoromethylbenzoyl)-cyclohexane-1,3-dione and 2-(2-chloro-4-methanesulfonylbenzoyl)-cyclohexane-1,3-dione. *Toxicol. Appl. Pharmacol.* **133(1)**, 12-19 (1995).
2. Grompe, M., Lindstedt, S., al-Dhalimy, M., *et al.* Pharmacological correction of neonatal lethal hepatic dysfunction in a murine model of hereditary tyrosinaemia type I. *Nat. Genet.* **10(4)**, 453-460 (1995).
3. Suzuki, Y., Oda, K., Yoshikawa, Y., *et al.* A novel therapeutic trial of homogentisic aciduria in a murine model of alkaptonuria. *J. Hum. Genet.* **44(2)**, 79-84 (1999).

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, 09/29/2022

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