

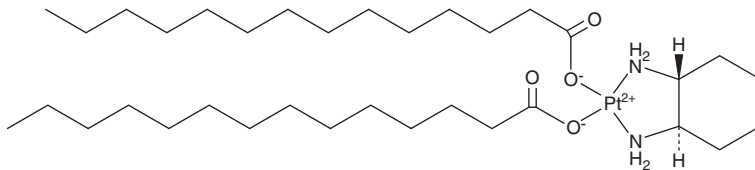
PRODUCT INFORMATION



Miriplatin

Item No. 43264

CAS Registry No.: 141977-79-9
Formal Name: (SP-4-2)-[(1R,2R)-1,2-cyclohexanediamine-κN¹,κN²]bis(tetradecanoato-κO)-platinum
Synonym: SM-11355
MF: C₃₄H₆₈N₂O₄Pt
FW: 764.0
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

Miriplatin is supplied as a solid. A stock solution may be made by dissolving the miriplatin in the solvent of choice, which should be purged with an inert gas. Miriplatin is slightly soluble (0.1-1 mg/ml) in chloroform.

Description

Miriplatin is a lipophilic derivative of cisplatin (Item No. 13119) containing myristic acid (Item No. 13351).¹ It decreases growth but increases serum levels of aspartate aminotransferase (AST) and alanine transaminase (ALT) in an AH109A-TC rat orthotopic hepatocellular carcinoma model when administered intra-arterially. Formulations containing miriplatin have been used in the treatment of hepatocellular carcinoma.

Reference

1. Kishimoto, S., Noguchi, T., Yamaoka, T., *et al.* Antitumor effects of a novel lipophilic platinum complex (SM-11355) against a slowly-growing rat hepatic tumor after intra-hepatic arterial administration. *Biol. Pharm. Bull.* **23**(3), 344-348 (2000).

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, 05/28/2025

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