

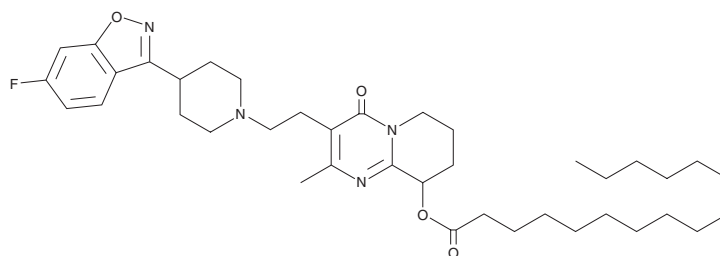
PRODUCT INFORMATION



Paliperidone Palmitate

Item No. 39644

CAS Registry No.: 199739-10-1
Formal Name: hexadecanoic acid, 3-[2-[4-(6-fluoro-1,2-benzisoxazol-3-yl)-1-piperidinyl]ethyl]-6,7,8,9-tetrahydro-2-methyl-4-oxo-4H-pyrido[1,2-a]pyrimidin-9-yl ester
Synonyms: 9-hydroxy Risperidone Palmitate, RO92670
MF: C₃₉H₅₇FN₄O₄
FW: 664.9
Purity: ≥98%
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

Paliperidone palmitate is supplied as a solid. A stock solution may be made by dissolving the paliperidone palmitate in the solvent of choice, which should be purged with an inert gas. Solvents such as methanol may be used. Paliperidone palmitate is slightly soluble in chloroform.

Paliperidone palmitate is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Paliperidone palmitate is a prodrug form of the atypical antipsychotic paliperidone (Item No. 15556).^{1,2} Formulations containing paliperidone palmitate have been used in the treatment of schizophrenia and schizoaffective disorder.

References

1. Chen, A., Cai, P., Peng, Y., *et al.* The role of alkyl chain length in the melt and solution crystallization of paliperidone aliphatic prodrugs. *IUCrJ* **11**(Pt. 1), 23-33 (2024).
2. Patel, H., Patel, P., Modi, N., *et al.* Rats and rabbits as pharmacokinetic screening tools for long acting intramuscular depots: Case study with paliperidone palmitate suspension. *Xenobiotica* **49**(4), 415-421 (2019).

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, 02/05/2024

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