

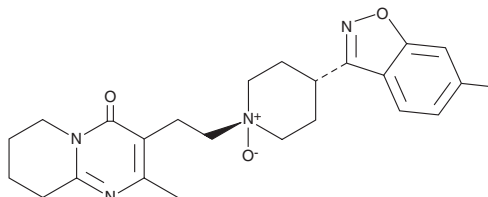
PRODUCT INFORMATION



Risperidone N-oxide

Item No. 46115

CAS Registry No.: 1894509-66-0
Formal Name: 3-[2-[*cis*-4-(6-fluoro-1,2-benzisoxazol-3-yl)-1-oxido-1-piperidinyl]ethyl]-6,7,8,9-tetrahydro-2-methyl-4H-pyrido[1,2-*a*]pyrimidin-4-one
Synonym: *cis*-Risperidone N-oxide
MF: C₂₃H₂₇FN₄O₃
FW: 426.5
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

Risperidone N-oxide is supplied as a solid. A stock solution may be made by dissolving the risperidone N-oxide in the solvent of choice, which should be purged with an inert gas. Risperidone N-oxide is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Risperidone N-oxide is a potential impurity found in commercial preparations of the atypical antipsychotic risperidone.^{1,2}

References

1. Tomar, R.S., Joseph, T.J., Murthy, A.S., *et al.* Identification and characterization of major degradation products of risperidone in bulk drug and pharmaceutical dosage forms. *J. Pharm. Biomed. Anal.* **36**(1), 231-235 (2004).
2. Germann, D., Kurylo, N., and Han, F. Risperidone. *Profiles Drug Subst. Excip. Relat. Methodol.* **37**, 313-361 (2012).

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, 08/11/2026

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