

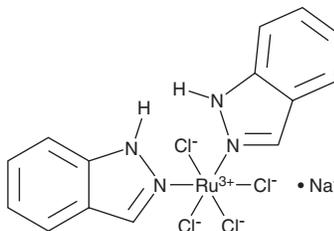
PRODUCT INFORMATION



NKP-1339

Item No. 35123

CAS Registry No.: 197723-00-5
Formal Name: (OC-6-11)-tetrachlorobis(1H-indazole- κ N²)-ruthenate(1-), monosodium salt
Synonym: KP1339
MF: C₁₄H₁₂Cl₄N₄Ru • Na
FW: 502.1
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

NKP-1339 is supplied as a solid. A stock solution may be made by dissolving the NKP-1339 in the solvent of choice, which should be purged with an inert gas. NKP-1339 is soluble in ethanol.

Description

NKP-1339 is a ruthenium complex with anticancer activity.¹ It is cytotoxic to HT-29 and SW480 cells (IC₅₀s = 24.46 and 123.3 μ M, respectively). NKP-1339 induces mitochondrial apoptosis in SW480 cells. *In vivo*, NKP-1339 (10.5, 21, and 42 mg/kg) reduces tumor growth in a Sal/N murine fibrosarcoma model.²

References

1. Kapitza, S., Pongratz, M., Jakupec, M.A., *et al.* Heterocyclic complexes of ruthenium(III) induce apoptosis in colorectal carcinoma cells. *J. Cancer Res. Clin. Oncol.* **131**(2), 101-110 (2005).
2. Hudej, R., Turel, I., Kanduser, M., *et al.* The influence of electroporation on cytotoxicity of anticancer ruthenium(III) complex KP1339 *in vitro* and *in vivo*. *Anticancer Res.* **30**(6), 2055-2063 (2010).

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, 12/09/2022

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