

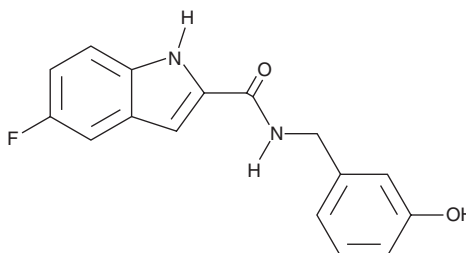
PRODUCT INFORMATION



KX1-004

Item No. 22198

CAS Registry No.: 518058-84-9
Formal Name: 5-fluoro-N-[(3-hydroxyphenyl)methyl]-1H-indole-2-carboxamide
MF: C₁₆H₁₃FN₂O₂
FW: 284.3
Purity: ≥98%
UV/Vis.: λ_{max}: 289 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

KX1-004 is supplied as a crystalline solid. A stock solution may be made by dissolving the KX1-004 in the solvent of choice. KX1-004 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of KX1-004 in these solvents is approximately 2.5, 20, and 30 mg/ml, respectively.

Description

KX1-004 is a non-ATP competitive inhibitor of Src protein tyrosine kinase (Src-PTK; IC₅₀ = 40 μM).¹ It protects against a permanent threshold shift in the auditory threshold and against outer hair cell loss in chinchillas following noise exposure at 106 dB when used at 30, 50, or 100 μM on the round window membrane of the ear. KX1-004 (50 mg/kg) is also effective against chronic noise exposure when administered systemically in chinchillas.²

References

1. Harris, K.C., Hu, B., Hangauer, D., *et al.* Prevention of noise-induced hearing loss with Src-PTK inhibitors. *Hear Res.* **208(1-2)**, 14-25 (2005).
2. Bielefeld, E.C., Hynes, S., Pryznosch, D., *et al.* A comparison of the protective effects of systemic administration of a pro-glutathione drug and a Src-PTK inhibitor against noise-induced hearing loss. *Noise Health* **7(29)**, 24-30 (2005).

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, 11/07/2022

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