

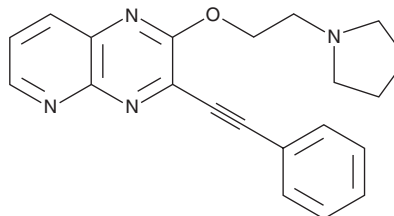
PRODUCT INFORMATION



GK921

Item No. 34359

CAS Registry No.: 1025015-40-0
Formal Name: 3-(2-phenylethynyl)-2-[2-(1-pyrrolidinyl)ethoxy]-pyrido[2,3-b]pyrazine
MF: C₂₁H₂₀N₄O
FW: 344.4
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 362, 378 nm
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

GK921 is supplied as a solid. A stock solution may be made by dissolving the GK921 in the solvent of choice, which should be purged with an inert gas. GK921 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of GK921 in these solvents is approximately 30 mg/ml.

GK921 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GK921 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. GK921 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GK921 is an inhibitor of transglutaminase 2 (TG2; IC₅₀ = 7.71 μM).¹ It is cytotoxic against a panel of eight renal cell carcinoma (RCC) cell lines (mean GI₅₀ = 0.905 μM). GK921 (8 mg/kg) suppresses tumor growth in ACHN and Caki-1 RCC mouse xenograft models.

Reference

1. Ku, B.M., Kim, S.-J., Kim, N., *et al.* Transglutaminase 2 inhibitor abrogates renal cell carcinoma in xenograft models. *J. Cancer Res. Clin. Oncol.* **140**(5), 757-767 (2014).

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, 11/22/2022

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