

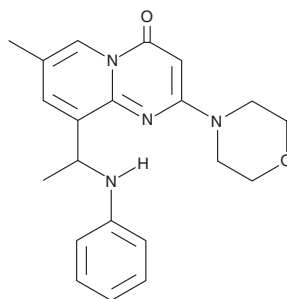
PRODUCT INFORMATION



TGX-221

Item No. 10007349

CAS Registry No.: 663619-89-4
Formal Name: 7-methyl-2-(4-morpholinyl)-9-[1-(phenylamino)ethyl]-4H-pyrido[1,2-a]pyrimidin-4-one
MF: C₂₁H₂₄N₄O₂
FW: 364.4
Purity: ≥98%
UV/Vis.: λ_{max}: 268 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

TGX-221 is supplied as crystalline solid. A stock solution may be made by dissolving the TGX-221 in the solvent of choice, which should be purged with an inert gas. TGX-221 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TGX-221 in ethanol is approximately 50 mg/ml and approximately 30 mg/ml in DMSO and DMF.

TGX-221 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TGX-221 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. TGX-221 has a solubility of approximately 0.5 mg/ml in a 1:8 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Phosphatidylinositol 3-kinase (PI3K) catalyzes the phosphorylation of phosphatidylinositol at the 3 position to produce the second messengers phosphatidylinositol-3,4-bisphosphate (PtdIns-(3,4)-P₂) and PtdIns-(3,4,5)-P₃.¹⁻³ Class 1 PI3Ks are composed of a p110 catalytic subunit, of which there are 4 isoforms (p110α, p110β, p110δ, and p110γ), and a p85 regulatory subunit.³ TGX-221 is a potent, selective, and cell permeable inhibitor of PI3K p110β.⁴ Inhibition appears to occur at the ATP binding site based on the observed increase in IC₅₀ value from 5 to ~50 nM at ATP concentrations of 50 μM and 1 mM, respectively. TGX-221 inhibits PtdIns-(3,4)-P₂ production in platelets with an IC₅₀ value of 50 nM.⁴ Selective inhibition of PI3K p110β results in defective platelet thrombus formation and defines PI3K as a target for antithrombotic therapy.⁴

References

1. Rameh, L.E. and Cantley, L.C. The role of phosphoinositide 3-kinase lipid products in cell function. *J. Biol. Chem.* **274**, 8347-8350 (1999).
2. Vivanco, I. and Sawyers, C.L. The phosphatidylinositol 3-kinase-AKT pathway in human cancer. *Nature Reviews Cancer* **2**, 489-501 (2002).
3. Hennessy, B.T., Smith, D.L., Ram, P.T., et al. Exploiting the PI3K/AKT pathway for cancer drug discovery. *Nature Reviews Drug Discovery* **4**, 988-1004 (2005).
4. Jackson, S.P., Schoenwaelder, S.M., Goncalves, I., et al. PI 3-kinase p110β: A new target for antithrombotic therapy. *Nature Med.* **11**(5), 507-514 (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, 01/03/2023

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