

PI-103

Catalog #: 27502-1

Lot #: 141111

Size: 5 mg

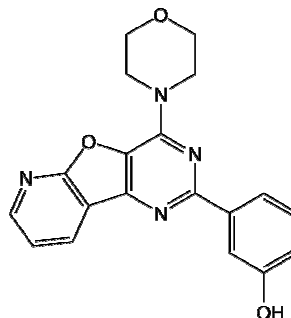
Structure:

CAS Registry #: 371935-74-9

Purity: ≥98%

Chemical Formula: C₁₉H₁₆N₄O₃

Molecular Weight: 348.36



Description: PI-103 is a multi-targeted inhibitor of PI3K for p110 α , p110 β , p110 γ , and p110 δ with reported IC₅₀ values of 2 nM, 3 nM, 3 nM, and 15 nM respectively. PI-103 is less potent towards mTOR/DNA-PK with reported IC₅₀ values of 30 nM and 23 nM respectively.

Appearance: Crystalline solid

Solubility: Soluble in DMSO or DMF at approximately 10 mg/ml. Poorly soluble in aqueous solutions. If necessary, first solubilize in DMSO or DMF and dilute into an aqueous solution, such as PBS.

Storage/Stability: At least 2 years as supplied at -20°C. At least 6 months in DMSO at -80°C. Do not store aqueous solutions for longer than one day. *Avoid freeze/thaw cycles.*

References:

1. Zou, Z.Q., *et al.*, *Int. J. Mol. Med.* 2009; **24(1)**: 97-101.
2. Lopez-Fauqued, M. *et al.*, *Int. J. Canc.* 2010; **126(7)**: 1549-1561.