

## PI-103

**Catalog #:** 27502-2

**Lot #:** 141111

**Size:** 25 mg

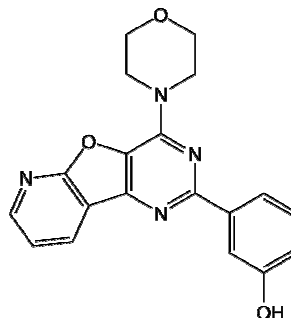
**Structure:**

**CAS Registry #:** 371935-74-9

**Purity:** ≥98%

**Chemical Formula:** C<sub>19</sub>H<sub>16</sub>N<sub>4</sub>O<sub>3</sub>

**Molecular Weight:** 348.36



**Description:** PI-103 is a multi-targeted inhibitor of PI3K for p110 $\alpha$ , p110 $\beta$ , p110 $\gamma$ , and p110 $\delta$  with reported IC<sub>50</sub> values of 2 nM, 3 nM, 3 nM, and 15 nM respectively. PI-103 is less potent towards mTOR/DNA-PK with reported IC<sub>50</sub> 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.