

## GDC-0941, Free Base

**Catalog #:** 27000

**Lot #:** 111107

**Size:** 10 mg

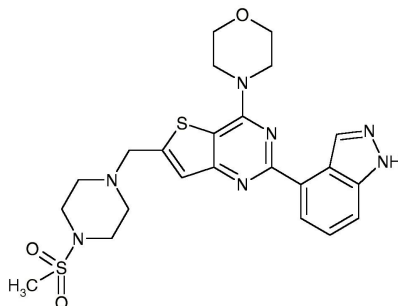
**Structure:**

**CAS Registry #:** 957054-30-7

**Purity:** ≥ 95%

**Chemical Formula:** C<sub>23</sub>H<sub>27</sub>N<sub>7</sub>O<sub>3</sub>S<sub>2</sub>

**Molecular Weight:** 513.64



**Description:** GDC-0941 is a phosphatidylinositol 3-kinase (PI3K) inhibitor.

**Appearance:** White solid.

**Solubility:** Soluble in DMSO at 66 mg/mL; very poorly soluble in ethanol; very poorly soluble in water; maximum solubility in plain water is estimated to be about 10-50 µM buffers, serum, or other additives may increase or decrease the aqueous solubility.

**Biological Activity:** GDC-0941 against p110α/p120γ IC<sub>50</sub>=0.003µM. The IC<sub>50</sub> for inhibiting cell proliferation in PC3 (prostate cancer) or U87MG (glioma) cells is 0.28 µM and 0.95 µM, respectively.

**Storage/Stability:** Store at or below -20 °C.

**Quality Control:** The purity was determined by HPLC.

### References:

1. Folkes, A.J. *et al. J. Med. Chem.* **51**: 5522-5532 (2008).
2. Yao, E. *et al Clin. Cancer Res.* **15**: 4147-4156 (2009).