

SGC707

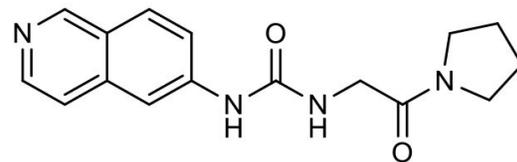
Catalog No: 14130, 14131**Format:** 25 mg, 5 mg**Chemical Properties:**

MW = 298.3

 $C_{16}H_{18}N_4O_2$

CAS 1687736-54-4

Physical Properties: Beige crystalline powder, MP = 233-235°C

Names: 1-Isoquinolin-6-yl-3-(2-oxo-2-pyrrolidin-1-yl-ethyl)-urea; SGC707

Structure of SGC707.

Pharmacology: Potent and selective allosteric inhibitor of protein arginine methyltransferase 3 (PRMT3)(ref 1). $IC_{50} = 31$ nM and selective against 31 other methyltransferases. An extremely useful probe for investigating the role of PRMT3 in normal and pathophysiology. Cell permeable and active *in vivo*.

Solubilization: May be dissolved in DMSO (35 mg/ml)**Fluorescent Properties:** N/A**Quality Control:**

>98% (TLC); NMR (Conforms)

References:

1. HU Kaniskan *et al. Angew.Chem. Int. Ed. Engl.* 2015, 54:5166

Storage and Guarantee: Store desiccated as supplied at room temperature for up to 2 years. Store solutions at -20°C for up to 3 months. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.