

Catalog No: 14096, 14097

Format: 50 mg, 10 mg

Chemical Properties:

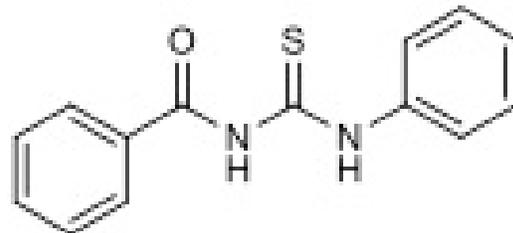
MW = 256.3

C₁₄H₁₂N₂OS

CAS 4921-82-8

Physical Properties: Beige powder

Names: 1-Benzoyl-3-phenyl-2-thiourea



Pharmacology: Highly potent and isozyme selective activator of recombinant HDAC8. Increases the activity of HDAC8 by a factor of 12 at 10 μM (ref1). No activation of other HDACs is observed. Rescues the activity of catalytically compromised HDAC8 mutants *in vitro* (ref2).

Solubilization: May be dissolved in DMSO (50 mg/ml); Ethanol (8 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (TLC); NMR (Conforms)

References:

1. RK Singh *et al. Bioorg. Med. Chem. Lett.* 2011, 21:5920
2. C Decroos *et al. ACS Chem. Biol.* 2014, 9:2157

Storage and Guarantee: Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.