

CPTH2

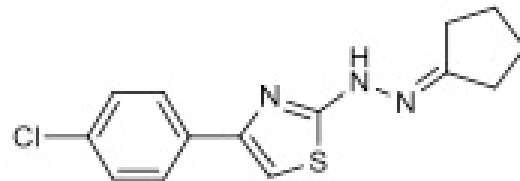
Catalog No: 14094, 14095**Format:** 25 mg, 5 mg**Chemical Properties:**

MW = 291.8

C₁₄H₁₄ClN₃S

CAS 357649-93-5

Physical Properties: White powder

Names: Cyclopentylidene-[4-(4'-chlorophenyl)thiazol-2-yl]hydrazone

Pharmacology: CPTH2 is a selective inhibitor of the histone acetyltransferase Gcn5P both *in vitro* and *in vivo* (IC₅₀ = 0.8 M) (ref1). *In vitro* the HAT inhibitory activity of CPTH2 was caused via bridge-interaction with histone H3. *In vivo* it decreased acetylation of bulk histone H3 at the H3K14ac site.

Solubilization: May be dissolved in DMSO (>25 mg/ml)**Fluorescent Properties:** N/A**Quality Control:**

>98% (TLC); NMR (Conforms)

References:

1. F Chimenti *et al. J. Med. Chem.* 2009, 52:530

Storage and Guarantee: Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month. 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.