

FK-866 HCl

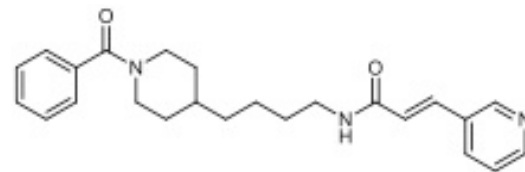
Catalog No: 14030, 14031**Format:** 25 mg, 5 mg**Chemical Properties:**

MW = 428.0

C₂₄H₂₉N₃O₂ HCl

CAS 658084-64-1

Physical Properties: Off white powder (hygroscopic), MP = 92-96°C

Names: FK-866 HCl

Chemical structure of FK-866.

Pharmacology: Nicotinamide phosphoribosyltransferase (Nampt) inhibitor (K_i = 0.4 nM). FK-866 is a highly specific non-competitive inhibitor of Nampt causing gradual NAD⁺ depletion (ref 1). Induces tumor cell apoptosis (ref 1) and causes premature senescence in normal human smooth muscle cells which may be linked to decreased activity of the NAD⁺-dependent enzyme SIRT1 (ref 2). Extends lifespan of human cells. Cell permeable.

Solubilization: May be dissolved in water (40 mg/ml)**Fluorescent Properties:** N/A**Quality Control:**

>98% (TLC); NMR (Conforms)

References:

1. M Hasmann *et al.* *Cancer Res.* 2003, 63:7436
2. E van der Veer *et al.* *J. Biol. Chem.* 2007, 282:10841

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