

ND-630

Catalog #: 27071-2

Lot #: 170509

Size: 5 mg

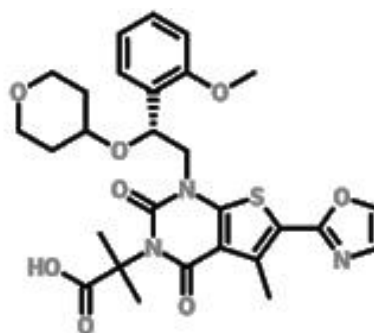
Structure:

CAS Registry #: 1434635-54-7

Purity: ≥99%

Chemical Formula: C₂₈H₃₁N₃O₈S

Molecular Weight: 569.63



Description: ND-630 (NDI 010976) is a highly specific, reversible inhibitor of ACC1 (IC₅₀ = 2 nM) and ACC2 (IC₅₀ = 6 nM). ND-630 interacts within the ACC phosphopeptide acceptor and dimerization site to prevent dimerization of ACC. Because the dimerization site is not conserved among other mammalian carboxylases, ND-630 lacks the ability to inhibit other mechanistically-related enzymes. In rats with diet-induced obesity, ND-630 reduced hepatic steatosis, improved insulin sensitivity, and reduced weight gain, suggesting ACC inhibition may be useful in treating metabolic disorders, including type 2 diabetes mellitus, and fatty liver disease.

Appearance: White powder

HPLC (Area %): 94.7

Storage/Stability: Store at -20°C for up to a year. Dissolve in anhydrous DMSO and store DMSO stock solutions at 4°C for a couple of days or store stock solutions at -20°C in aliquots for up to a month. Avoid freeze/thaw cycles of solutions.

References:

1. Harriman, G., *et al.*, *PNAS*. 2016; **113 (13)**: E1796-E1805.