

Danoprevir (ITMN-191)

Catalog #: 27052

Lot #: 130909

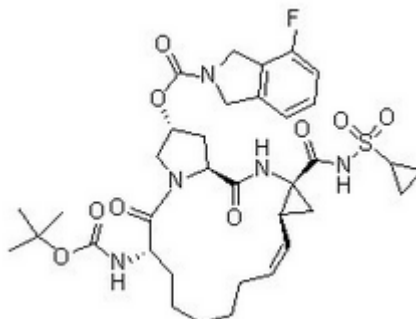
Size: 5 mg

Structure:

CAS Registry #: 850876-88-9

Purity: ≥ 95%

Chemical Formula: C₃₅H₄₆FN₅O₉S



Molecular Weight: 731.83

Description: Danoprevir is a peptidomimetic inhibitor. It inhibits of the NS3/4A protease of hepatitis C virus (HCV).

Appearance: Off-white powder.

Solubility: Soluble in DMSO at 144 mg/mL, in ethanol at 144 mg/mL, and poorly soluble in H₂O.

Biological Activity: Danoprevir inhibits NS3/4A protease of HCV with an IC₅₀ of 0.2-3.5 nM. It also shows antiviral inhibition effects against HCV genotypes 1, 4 and 6 with IC₅₀ of 2-3 nM.

Storage/Stability: Room temperature for 24 months as powder, +4°C for 2 weeks in DMSO, or -80°C for 6 months in anhydrous DMSO. Keep DMSO stock solutions out of contact with air. Protect from moisture.

Quality Control: The purity was determined by HPLC.

References:

1. Robinson M, *et al.*, *Curr Protoc Microbiol*; 17,17.7 (2011).
2. Saeed M., *et al.*, *J Antimicrob Agents Chemother*; **56(10)**: 5365-5373 (2012).