

VX-222 (VCH-222)

Catalog #: 27053

Lot #: 130909

Size: 5 mg

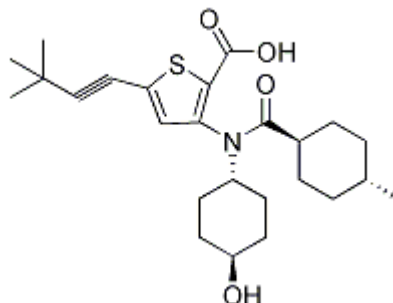
Structure:

CAS Registry #: 1026785-55-6

Purity: ≥ 99%

Chemical Formula: C₂₅H₃₅NO₄S

Molecular Weight: 445.61



Description: VX-222 (VCH-222) is a potent, selective, novel inhibitor of non-nucleoside polymerase, specifically the HCV RNA-dependent RNA. It works by inhibiting HCV NS5B of genotype 1a and 1b, the replication of subgenomic HCV genotype 1a and 1b, the 1b/Con1 HCV subgenomic replicon.

Appearance: White to off-white solid or powder.

Solubility: Soluble in DMSO at 89 mg/mL, in ethanol at 89 mg/mL, and poorly soluble in H₂O.

Biological Activity: : VX-222 (VCH-222) is shown Inhibit HCV NS5B of genotype 1a and 1b (with IC₅₀ of 0.94 and 1.2 μM, respectively), the replication of subgenomic HCV genotype 1a and 1b (with an EC₅₀ of 22.3 and 11.2 nM, respectively.), and the 1b/Con1 HCV subgenomic replicon at EC₅₀ of 5 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. Bedard J, *et al.*, *J.Hepatol*; **50**: S340 (2009).
2. Chouret N., *et al.*, *J.Hepatol*; **50**: S341-S342 (2009).
3. Yi, G., *et al.*, *Antimicrob Agents Chemother*; **56(2)**: 830 – 837 (2012).