

BMS-790052

Catalog #: 27054

Lot #: 130906

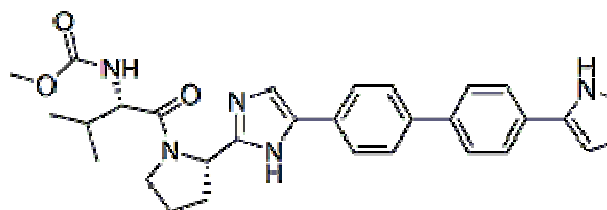
Size: 5 mg

Structure:

CAS Registry #: 1009119-64-5

Purity: ≥ 97%

Chemical Formula: C₄₀H₅₀N₈O₆



Molecular Weight: 738.88

Description: BMS-790052 is a highly potent, selective inhibitor of HCV NS5A. It allows for inhibition of a broad range of HCV replicon genotypes as well as the JFH-1 genotype 2a infectious virus in cell culture.

Appearance: A white crystalline powder.

Solubility: Soluble in DMSO at 148 mg/ml, in ethanol at 148 mg/ml, and poorly soluble in H₂O.

Biological Activity: BMS-790052 is a highly potent, selective inhibitor of HCV with an EC₅₀ of 9 and 50 pM for the HCV genotypes 1b and 1a, respectively.

Storage/Stability: Room temperature for 24 months as powder, +4°C for 2 weeks in DMSO, or -70°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. Bush CO., *et al.*, *PLoS One*. **8(6)**:e65273 (2013).
2. Kumthip K., *et al.*, *J Virol*. **86(16)**:8581-8591(2012).