

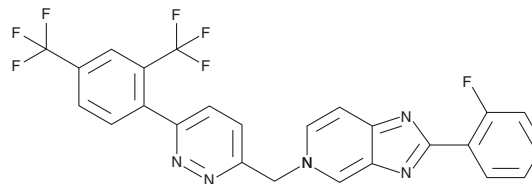
# PRODUCT INFORMATION



## Tegobuvir

Item No. 36603

**CAS Registry No.:** 1000787-75-6  
**Formal Name:** 5-[[6-[2,4-bis(trifluoromethyl)phenyl]-3-pyridazinyl]methyl]-2-(2-fluorophenyl)-5H-imidazo[4,5-c]pyridine  
**Synonyms:** GS-9190, GS-333126  
**MF:** C<sub>25</sub>H<sub>14</sub>F<sub>7</sub>N<sub>5</sub>  
**FW:** 517.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 243, 301 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Tegobuvir is supplied as a solid. A stock solution may be made by dissolving the tegobuvir in the solvent of choice, which should be purged with an inert gas. Tegobuvir is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

### Description

Tegobuvir is a non-nucleoside inhibitor of hepatitis C virus (HCV) non-structural protein 5B (NS5B).<sup>1</sup> It reduces viral infectivity in cells expressing wild-type HCV 1b NS5B but not mutant NS5Bs. Tegobuvir selectively inhibits replication of HCV genotypes 1a and 1b over genotype 2a in Huh7-Lunet cells (EC<sub>50</sub>s = 13.8, 0.8, and 21,900 nM, respectively).

### Reference

1. Shih, I.-H., Vliegen, I., Peng, B., *et al.* Mechanistic characterization of GS-9190 (Tegobuvir), a novel nonnucleoside inhibitor of hepatitis C virus NS5B polymerase. *Antimicrob. Agents Chemother.* **55**(9), 4196-4203 (2011).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/20/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)