

Vismodegib (GDC-0449)

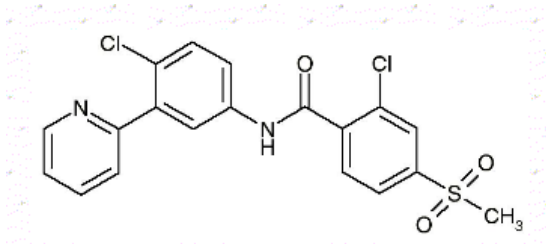
Catalog #: 27015

Lot #: 170124

Size: 50 mg

Structure:

CAS Registry #: 879085-55-9



Purity: ≥ 99%

Chemical Formula: C₁₉H₁₄Cl₂N₂O₃S

Molecular Weight: 421.30

Description: Vismodegib (GDC-0449) is a potent hedgehog (Hh) signaling pathway inhibitor.

Appearance: A crystalline solid.

Solubility: Soluble in DMSO at 200 mg/mL; soluble in ethanol at 10 mg/mL with warming; very poorly soluble in water; maximum solubility in plain water is estimated to be about 10-20 μM; buffers, serum, or other additives may increase or decrease the aqueous solubility.

Biological Activity: Inhibits two ABC transporters, ABCG2/BCRP and ABCB1/Pgp, with IC₅₀ values of 1.4 and 3.0 μM, respectively.

Storage/Stability: Store at or below -20°C.

Quality Control: The purity was determined by HPLC.

References:

1. Von Hoff, D.D., *et al.* *N. Engl. J. Med.* **361**: 1164-1172 (2009).
2. Zhang, Y., *et al.* *Neoplasia* **11**: 96-101 (2009).