

Product Information

Description

ICG 1 is a small molecule inhibitor that targets Wnt/ β -catenin pathway. It inhibits TCF/ β -catenin mediated transcription by competing with β -catenin for CBP (CREB binding protein) but not p300 binding with an IC₅₀ of 3 μ M. It can be used to specifically evaluate the role of CBP/ β -catenin association on WNT signaling and cell physiology.

Target:

CBP

Chemical Name:

(6S,9aS)-N-benzyl-6-[(4-hydroxyphenyl)methyl]-8-(naphthalen-1-ylmethyl)-4,7-dioxo-3,6,9,9a-tetrahydro-2H-pyrazino[1,2-a]pyrimidine-1-carboxamide

Molecular Formula:C₃₃H₃₂N₄O₄**Supplied As:**

Solid

Solubility:

≥27.43 mg/mL in DMSO; insoluble in H₂O; ≥35.47 mg/mL in EtOH with ultrasonic. For obtaining a higher solubility, please warm the tube at 37°C and shake it in the ultrasonic bath for a while.

Stability:

At least 6 months at -20°C. Avoid freeze/thaw cycles.

Storage:

-20°C

MW:

548.63 Da

Cas No.

847591-62-2

Purity:

≥98%

Chemical Structure

