

## Product Information

## Description

ICG 1 is a small molecule inhibitor that targets Wnt/ $\beta$ -catenin pathway. It inhibits TCF/ $\beta$ -catenin mediated transcription by competing with  $\beta$ -catenin for CBP (CREB binding protein) but not p300 binding with an IC<sub>50</sub> of 3  $\mu$ M. It can be used to specifically evaluate the role of CBP/ $\beta$ -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:

C<sub>33</sub>H<sub>32</sub>N<sub>4</sub>O<sub>4</sub>

## Supplied As:

Solid

## Solubility:

$\geq 27.43$  mg/mL in DMSO; insoluble in H<sub>2</sub>O;  $\geq 35.47$  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:

 $\geq 98\%$ 

## Chemical Structure

